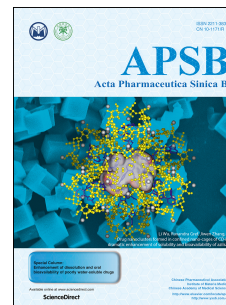


Journal Pre-proof

Inhibition of intracellular proton-sensitive Ca^{2+} -permeable TRPV3 channels protects against ischemic brain injury

Xiaoling Chen, Jingliang Zhang, KeWei Wang



PII: S2211-3835(22)00013-2

DOI: <https://doi.org/10.1016/j.apsb.2022.01.001>

Reference: APSB 1304

To appear in: *Acta Pharmaceutica Sinica B*

Received Date: 13 May 2021

Revised Date: 18 November 2021

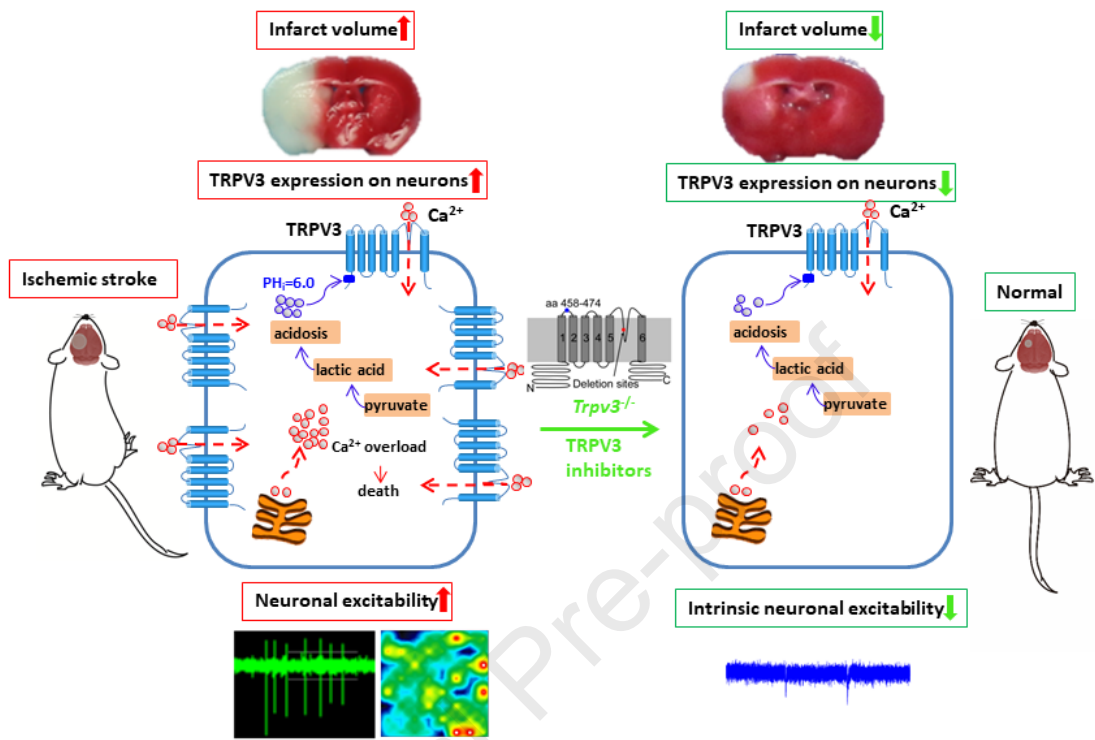
Accepted Date: 4 January 2022

Please cite this article as: Chen X, Zhang J, Wang K, Inhibition of intracellular proton-sensitive Ca^{2+} -permeable TRPV3 channels protects against ischemic brain injury, *Acta Pharmaceutica Sinica B*, <https://doi.org/10.1016/j.apsb.2022.01.001>.

This is a PDF file of an article that has undergone enhancements after acceptance, such as the addition of a cover page and metadata, and formatting for readability, but it is not yet the definitive version of record. This version will undergo additional copyediting, typesetting and review before it is published in its final form, but we are providing this version to give early visibility of the article. Please note that, during the production process, errors may be discovered which could affect the content, and all legal disclaimers that apply to the journal pertain.

© 2022 Chinese Pharmaceutical Association and Institute of Materia Medica, Chinese Academy of Medical Sciences. Production and hosting by Elsevier B.V. All rights reserved.

Graphical abstract



Original article

Inhibition of intracellular proton-sensitive Ca²⁺-permeable TRPV3 channels protects against ischemic brain injuryXiaoling Chen^{a,†}, Jingliang Zhang^{a,†}, KeWei Wang^{b,*}^a*Department of Molecular and Cellular Pharmacology, State Key Laboratory of Natural and Biomimetic Drugs, School of Pharmaceutical Science, Peking University, Beijing 100191, China*^b*Department of Pharmacology, Qingdao University School of Pharmacy, Qingdao 266021, China*

Received 26 August 2021; received in revised form 18 November 2021; accepted 2 December 2021

*Corresponding author. Tel.: +86 532 82991070.

E-mail address: wangkw@qdu.edu.cn (KeWei Wang).

†These authors made equal contributions to this work.

Running title: A causative role of overactive TRPV3 in brain stroke

Abstract Ischemic brain stroke is pathologically characterized by tissue acidosis, sustained calcium entry and progressive cell death. Previous studies focusing on antagonizing *N*-methyl-D-aspartate (NMDA) receptors have failed to translate any clinical benefits, suggesting a non-NMDA mechanism involved in the sustained injury after stroke. Here, we report that inhibition of intracellular proton-sensitive Ca²⁺-permeable transient receptor potential vanilloid 3 (TRPV3) channels protects against cerebral ischemia/reperfusion (I/R) injury. TRPV3 expression is upregulated in mice subjected to cerebral I/R injury. Silencing of TRPV3 reduces intrinsic neuronal excitability, excitatory synaptic transmissions, and also attenuates cerebral I/R injury in mouse model of transient middle cerebral artery occlusion (tMCAO). Conversely, overexpressing or re-expressing TRPV3 increases neuronal excitability, excitatory synaptic transmissions and aggravates cerebral I/R injury. Furthermore, specific inhibition of TRPV3 by natural forsythoside B decreases neural excitability and attenuates cerebral I/R injury. Taken together, our findings for the first time reveal a causative role of neuronal TRPV3 channel in progressive cell death after stroke, and blocking overactive TRPV3 channel may provide therapeutic potential for ischemic brain injury.

KEY WORDS TRPV3; Ca²⁺ influx; Acidosis; Cerebral ischemia/reperfusion injury; Excitotoxicity; Neural excitability; Forsythoside B; tMCAO; OGD; TRP

1. Introduction

As a result from an interruption of blood supply, ischemic brain injury is characterized by histopathological acidosis of brain tissues^{1,2}. Oxygen and glucose deprivation necessitates a switch

from aerobic metabolism to anaerobic glycolysis, which leads to the generation of lactic acid and intracellular protons^{3,4}. Accumulation of lactic acid as a byproduct of glycolysis and protons produced by adenosine triphosphate (ATP) hydrolysis causes a typical fall of intracellular pH between 6.5 to 6.0 in the ischemic area of the brain⁵⁻⁷. Consequently, tissue acidosis aggravates ischemic brain injury partly through acid-sensing ion channels⁸⁻¹⁰ and Na⁺/H⁺ exchanger isoform 1^{11,12}.

Despite the progress made in prevention, treatment for ischemic brain stroke still remains a major global challenge. The NMDA-based neuronal excitotoxicity theory is based on the observations that occurrence of cerebral ischemic injury causes the release of neuronal excitatory neurotransmitter glutamate that activates ionotropic glutamate receptors (iGluRs) and leads to an overload of Ca²⁺ influx, ultimately further aggravating ischemic injury and neuronal apoptosis and necrosis¹³⁻¹⁶. Over the past three decades, all NMDA receptor antagonists, however, have been unable to translate any clinical benefits for stroke treatment due to poor clinical outcomes or severe adverse reactions¹⁷⁻²⁰, thus suggesting that glutamate-independent mechanisms for calcium overload during ischemia may be involved in the neuronal death after stroke²¹⁻²³.

We have previously demonstrated that the warm-temperature sensitive and Ca²⁺-permeable cation transient receptor potential vanilloid 3 (TRPV3) channel is activated by intracellular protons with half-maximal activation (pH_{0.5}) at pH 6.1²⁴. The proton-mediated activation of TRPV3 causes an overload of Ca²⁺ influx, leading to cell death²⁴. TRPV3 is highly expressed in the brain and spinal cord, dorsal root ganglia (DRG)²⁵⁻²⁷. We therefore hypothesized that brain tissue acidosis during cerebral ischemia may trigger excessive activation of TRPV3 channels through intracellular rise of protons, which further causes intracellular Ca²⁺ overload and progressive cell death after ischemic injury.

To test this hypothesis, we investigated the role of intracellular proton-activated Ca²⁺-permeable TRPV3 channels in the pathogenesis of cerebral ischemia/reperfusion (I/R) injury. We observed an upregulation of TRPV3 expression in mouse brain subjected to I/R injury. Overexpressing or re-expressing TRPV3 increases neuronal firing and aggravates cerebral I/R injury. Silencing or pharmacological inhibition of TRPV3 suppresses neural excitability and also protects against cerebral I/R injury. Our findings reveal a non-NMDA-mediated neural excitotoxicity involving overactive TRPV3 in the progression of ischemic injury, and that blocking overactive proton-sensitive and Ca²⁺-permeable TRPV3 channel may provide an effective recovery and potential therapy for ischemic brain stroke.

2. Materials and methods

2.1. Chemicals

Forsythoside B (FB, molecular weight 756.7, 99% purity) was purchased from Shanghai Tongtong biotechnology Co., Ltd. (Shanghai, China). Antioxidant edaravone or redicava (Simcere) was used as a positive control for ischemia experiments. *N*-Methyl-D-aspartate (NMDA) receptor antagonist D-2-amino-5-phosphonovalerate (AP5, 50 $\mu\text{mol/L}$), α -amino-3-hydroxyl-5-methyl-4-isoxazole-propionate (AMPA)/kainate glutamate receptor antagonist 6-cyano-7-nitroquinoxaline-2,3-dione (CNQX, 10 $\mu\text{mol/L}$) were purchased from Sigma–Aldrich. Neurobiotin tracer *N*-(2-aminoethyl) biotinamide hydrochloride (Vector Laboratories), Alexa Fluor 488 or 647-conjugated streptavidin (Molecular Probes), sodium channel blocker tetrodotoxin (TTX, 500 nmol/L, Baomanbio), γ -aminobutyric acid A (GABA_A) receptor antagonist bicuculline (10 $\mu\text{mol/L}$, Abcam), and GABA_A receptor agonist muscimol (1 $\mu\text{mol/L}$, Abcam) were prepared before use, and the purity of each standard compound is no less than 98% by high-performance liquid chromatography analysis. FB and edaravone used in behavior experiments were diluted in sterile normal saline for intravenous injection.

2.2. Animals

Male adult C57BL/6J mice (10 ± 2 weeks) and KM mice (10 ± 2 weeks) of SPF grade were purchased from Charles River Laboratories (Beijing, China; permit No. SCXK 2012-0001). All animals were kept on a 12 h:12 h light–dark cycle with free access to food and water. Animals were handled in strict accordance with the “Guide for the Care and Use of Laboratory Animals” and the “Principles for the Utilization and Care of Vertebrate Animals”. All animal protocols were approved by the Institutional Animal Care and Use Committee at Peking University (Beijing, China) or Qingdao University (Qingdao, China) and efforts were made to minimize animal suffering and the number of animals used. *In vivo* experiments were blind to viral or drug treatment conditions during transient middle cerebral artery occlusion (tMCAO) and behavioral testing. C57BL/6J mice were used to generate the model of tMCAO, and KM mice were used to evaluate the neuroprotection of FB on ischemia injury.

Trpv3 deficient mice (Allele Symbol: *Trpv3*^{tm1Apat}, 10 ± 2 weeks) from kind gifts of Dr. Yong Yang were originally produced in the laboratory of Dr. Ardem Patapoutian²⁸ and bred in our animal facility. The mouse genotype was determined using PCR analysis. Briefly, genomic DNA was extracted from 0.2 cm tail snips using the alkali extraction method²⁹. A PCR reaction was then performed using Ex Taq polymerase (TaKaRa) and the following primers:

primer 1, 5'-GTAGAGTCAGGGCCCTCAGAGGAGCC-3';

primer 2, 5'-CTAATGTTGCAGGTACTGTGTCGCCCC-3';

primer 3, 5'-GTGGGCTCTATGGCTTCTGAGGCGG-3'.

The PCR reaction consisted of an initial 2 min at 95 °C, followed by 40 cycles of 30 s at 95 °C, 30 s at 59 °C, and 40 s at 72 °C. After the last cycle, the reaction was kept at 72 °C for 5 min and then held at 4 °C. A 304 bp band was observed for mice containing the wild-type (WT) allele, whereas a 182 bp band was seen for mice containing the mutant allele and heterozygotes contained both alleles. Male homozygotes *Trpv3*^{-/-} mice (10 ± 2 weeks) and WT were used to generate ischemia injury.

2.3. Generation of cerebral ischemia/reperfusion (I/R) injury by transient middle cerebral artery occlusion (tMCAO) in mice

Mice were anesthetized by intraperitoneal injection of pentobarbital sodium (60 mg/kg). After sterilizing with 75% alcohol, mouse scalp was cut and exposed to the skull. A probe was connected to the left skull by biomedical glue to monitor relative Local cerebral blood flow (LCBF) with laser Doppler measurement (Perimed). Transient focal cerebral ischemia was induced by tMCAO surgery³⁰. Briefly, a 6-0 nylon monofilament suture blunted at the tip and coated with silicone was introduced ~0.8 mm into the internal carotid artery *via* the external carotid artery, to occlude the origin of the left middle cerebral artery (MCA) in the circle of Willis. The LCBF should drop below 80% of the baseline at ischemia time during the surgery³¹. Reperfusion *via* MCA was allowed after 1.5 h by monofilament removal and the LCBF should recover over 80% of the baseline after reperfusion. Body temperature was maintained and monitored at 37 ± 0.5 °C by a heating operation table during the surgery and the reperfusion. For sham operations, the filament is inserted to occlude MCA before withdrawn immediately to allow instant reperfusion. The subsequent operation is identical to animals undergoing cerebral ischemia.

2.4. Neurologic deficit scoring

Neurologic deficit scores were evaluated at 24 h of reperfusion according to Longa's five-point scale method³². In detail, 0, no neurologic deficit; 1, a mild focal neurologic deficit: failure to extend left forepaw fully; 2, a moderate focal neurologic deficit: circling to the left; 3, a severe focal deficit: falling to the left; 4, mice did not walk spontaneously and had a depressed level of consciousness.

An expanded seven-point scale is used for evaluation of transgenic mice³⁰ on Days 1, 3 and 5. That is, 0: normal; 1: failure to extend left forepaw fully; 2: mild circling behavior with or without inconsistent rotation when picked up by the tail, <50% attempts to rotate to the contralateral side; 3:

mild consistent circling, >50% attempts to rotate to the contralateral side; 4: consistent strong and immediate circling, the mouse holds a rotation position for more than 1–2 s, with its nose almost reaching its tail; 5: severe rotation progressing into barreling, loss of walking or righting reflex; 6: comatose or moribund. The observer was blind to animal treatment, and one mouse obtaining no less than 2 scores is counted as a valid model.

2.5. Infarct volume measurement

Infarct volume was determined 24 h after reperfusion. Mice were perfused with saline to remove blood in brain tissue under anesthetizing. The brains were quickly removed, sectioned coronally into 6 slices at ~1-mm thickness and stained by immersion in the vital dye 1% 2,3,5-triphenyltetrazolium hydrochloride (TTC, Amresco, w/v) in phosphate-buffered saline (PBS) at 37 °C each side for 10 min^{33,34}, followed by immersion-fixation in 4% paraformaldehyde. Then TTC-stained sections were photographed and the extents of the normal and infarcted areas were analyzed using Adobe Photoshop CC and determined by the indirect method, which corrects for edema^{33,35}. The percentage of the corrected infarct volume was calculated by dividing the infarct volume by the total contralateral hemispheric volume to compensate for the effect of brain edema³⁶.

2.6. Measurement of brain water content

Brain water content was determined according to the wet-dry methods³⁷. In brief, brains were immediately weighed to obtain wet weight (W_1), then dried in an oven at 120 °C for 8 h and weighed again to obtain the dry weight (W_2). Brain water content was calculated as Eq. (1):

$$\text{Brain water content (\%)} = (W_1 - W_2) / W_1 \times 100 \quad (1)$$

2.7. Oxygen–glucose deprivation (OGD) in primary culture of mouse cortical neurons

Cortical neurons were removed from newborn mice (<24 h) and dissociated by 0.25% trypsinization³⁸. Cells were suspended in high-glucose DMEM (Gibco) containing 10% fetal bovine serum (Gibco), and plated on poly-D-lysine hydrobromide (Sigma–Aldrich) coated culture plates or 15 mm confocal culture dish at a density of about 1.2×10^5 cells/cm² for different experiments. After 4–6 h seeding, the medium was changed to phenol red-free neurobasal A medium (Gibco) supplemented with 2% B27 (Gibco), containing 0.5 mmol/L GlutaMAX- I, 0.5% penicillin–streptomycin, then one-half medium was refreshed every three days. At the 3rd day, to suppress the proliferation of glial cells, cytarabine (2.5 µg/mL) was added. Cells were maintained at 37 °C in a humidified atmosphere containing 95% air and 5% CO₂ and used for experiments after 7–9

days *in vitro*.

Seven to nine days *in vitro* neurons were prepared and the OGD model was carried out as described previously³⁹. The original media were removed, the cells were washed with 37 °C glucose-free and phenol red-free DMEM twice, and placed with a glucose-free and phenol red-free DMEM containing 10 mmol/L of sodium dithionite (Na₂S₂O₄), a deoxygenated reagent. A duration of 30 or 20 min later, the media were removed, the cells were washed with 37 °C neurobasal A medium twice, and returned to their original culture condition and maintained for 24 h until the assay of cell injury. FB (0.1, 1, and 10 μmol/L) and edaravone (10 μmol/L) were added to culture medium 30 min before exposed to OGD and lasted for reperfusion 24 h.

2.8. Cell viability and lactate dehydrogenase (LDH) assays

A cell counting kit-8 assay (CCK-8, Dojindo) was used to estimate cell viability. Briefly, a highly water-soluble tetrazolium salt, WST-8, is reduced by dehydrogenase activities in cells to give a yellow-color formazan dye, which is soluble in the tissue culture media. The amount of the formazan dye, generated by the activity of dehydrogenases in cells, is directly proportional to the number of living cells. CCK-8 reagents (10%) were added into wells, and the absorbance was measured at 450 nm using a microplate reader at 2 h after CCK-8 addition. At least three wells were measured for cell viability in each group⁴⁰.

The release of lactate dehydrogenase (LDH) into the culture medium was measured using LDH assay reagent (Promega) following the instructions of the kit menu. Briefly, the supernatant medium (50 μL) was transferred to 96-well plates, mixed with 50 μL of reaction solution in kits and incubated for 30 min at room temperature. Stop solution in kits was then added and the optical density was measured at 490 nm.

2.9. Stereotactic injection and histological examinations

Viral AAV2 constructs of mTRPV3 with enhanced green fluorescent protein (mTRPV3-EGFP) and control-EGFP carrying human synapsin (hsyn) promoter for neuron-specific expression were designed and packed by Shanghai Hanbio Biotechnology Co., Ltd. (Shanghai, China). For viral injection, mice were anesthetized with pentobarbital sodium (60 mg/kg, intraperitoneal injection, i.p.) and placed in a stereotaxic apparatus (RWD). Mice were bilaterally injected with AAV virus (~10¹² infections units per mL) into the cortex (coordinates from bregma: +0.5 mm anterior/posterior, ±2.0 mm medial/lateral, -1.8 mm dorsal/ventral, 2 μL per point) and the striatum (coordinates from bregma: +0.5 mm anterior/posterior, ±2.0 mm medial/lateral, -3.5 mm dorsal/ventral, 2 μL per point)

at a slow rate (~100–150 nL/min). The injection needle was slowly withdrawn 5 min after the virus infusion. For the left tMCAO surgery, viruses were injected in the left cortex and striatum. tMCAO experiments or electrophysiological recordings were performed at least 30 days after viral injection.

After 30 days of viral injection, mice were deeply anesthetized and perfused with normal saline followed by 4% paraformaldehyde in PBS. Brains were post-fixed overnight in the same 4% paraformaldehyde solution and dehydrated by immersion in 20% and 30% sucrose solution in PBS for 1 day. Frozen brains were sectioned at 20 µm with a sliding microtome (Leica) in the coronal plane. Brain slices from mice injected with AAV viruses were directly examined under a fluorescent microscope.

2.10. Immunostaining and confocal microscopy

HEK 293T cells were transfected with mouse-TRPV3-EGFP plasmid using Lipofectamine 2000 (Invitrogen) following the manufacturer's instructions and were maintained at 37 °C under 5% CO₂ in DMEM supplemented with 10% fetal bovine serum. HEK 293T cells expressing TRPV3 and mouse cortical neurons in 15 mm confocal culture dish were fixed in 4% paraformaldehyde for 15 min at 4 °C after removal of culture medium. The cells were added with 0.3% Triton X-100 (Amresco) in PBS for 15 min and blocked by 10% sheep serum in 0.01 mol/L PBS for 1 h at room temperature before incubated with primary antibodies overnight at 4°C. The rabbit monoclonal NeuN antibody (1:1000; Abcam), mouse monoclonal TRPV3 antibody (1:1000; Abcam). The immunoreactivity was visualized with Alexa Fluor 488- or 594-conjugated secondary antibodies (1:200; ZSGB-BIO). Immunocytochemistry was used for visualization of TRPV3 proteins in cells using a confocal microscope (Leica Microsystems).

2.11. RNA isolation, reverse transcription, qPCR analysis

After MCAO surgery at 6, 12, and 24 h, mouse brains were divided into ischemia part and contralateral part for extraction of RNAs. Total RNAs were extracted with TRIzol reagent (Sigma) from mouse cerebral tissues according to the manufacturer's instructions. RNA (4 µg) was subjected to reverse transcription using a GoScript Reverse Transcription System (Promega), and resulting cDNAs were subjected to quantitative PCR analysis using the GoTaq qPCR master mix (Promega) and specific primers in a 7500 Fast Real-Time PCR System (Applied Biosystems). PCR primer sequences were as follows:

mouse *Trpv3* forward 5'-ACAGGTTTCATCAACGCTGAGT-3';

mouse *Trpv3* reverse 5'-CACCCGCTGCTATAAGCACT-3';

mouse *Actb* forward 5'-GGCTGTATTCCCCTCCATCG-3';

mouse *Actb* reverse 5'-CCAGTTGGTAACAATGCCATGT-3';

mGapdh-forward 5'-GTCTTCACCACCATGGAGA-3';

mGapdh-reverse 5'-AAGCAGTTGGTGGTGCAG-3'.

Gapdh levels, *Actb* levels and *Trpv3* mRNA expression in contralateral cortex were used as an endogenous control for normalization using the $\Delta\Delta Ct$ method⁴¹. In brief, test (T): $\Delta Ct^T = [Ct^T \text{ (target gene)} - Ct^T \text{ (internal control)}] - [Ct^T \text{ (contralateral target gene)} - Ct^T \text{ (contralateral internal control)}]$; Sham control (C): $\Delta Ct^C = [Ct^C \text{ (target gene)} - Ct^C \text{ (internal control)}] - [Ct^C \text{ (contralateral target gene)} - Ct^C \text{ (contralateral internal control)}]$; $\Delta\Delta Ct = \Delta Ct^T - \Delta Ct^C$; Amount of the target = $2^{-\Delta\Delta Ct}$.

2.12. Western blot

Brain regions from the right and left hemispheres corresponding to the ischemic core, penumbra and contralateral cerebral were dissected^{42,43}. Briefly, a midline between the two hemispheres was identified before a longitudinal cut (from top to bottom) approximately 2 mm from the midline through each hemisphere. A transverse diagonal cut was made at approximately “10 o’clock” position to separate the penumbra (adjacent cortex) from the core (*i.e.*, striatum and overlying cortex). We made minor modifications for estimates of core and penumbra based on our earlier investigations by TTC staining. Proteins were extracted from cerebral regions in cold RIPA Lysis Buffer containing 2% Cocktail (Roche). Protein samples were loaded on 8% sodium dodecyl sulfate-polyacrylamide gel electrophoresis before transferred using gel electrophoresis to polyvinylidene fluoride (PVDF) membranes (Millipore). After 5% milk blocking, PVDF membranes were incubated with mouse monoclonal TRPV3 antibody or mouse β -actin antibody (1:5000, Abcam) at 4 °C overnight. The membranes were then incubated with their corresponding secondary horseradish peroxidase-conjugated antibodies and detected using an ECL Western blotting detection system (Millipore). The immunoreactive bands were scanned by Tanon 5200 instrument (Tanon), captured by the Tanon MP system (Tanon) and quantitatively analyzed by densitometry with Tanon GIS software (Tanon) or ImageJ software. For increase of low TRPV3 protein signal in Western blot assay, a reagent called “SuperSignal Western blot enhancer (Thermo Fisher)” was used. We used TRPV3 proteins as positive controls (TRPV3 protein expression on HEK293 cell line or on mouse skin) and checked molecular weight 91 kDa to make sure the TRPV3 protein band site.

2.13. Acute brain slice preparations and electrophysiological whole-cell recordings

Mice were anesthetized by intraperitoneal injection of pentobarbital sodium (60 mg/kg) and

decapitated before brains were dissected into ice-cold slice solution containing the following (in mmol/L): 110 choline chloride, 2.5 KCl, 1.25 NaH₂PO₄, 25 NaHCO₃, 0.5 CaCl₂, 7 MgCl₂, 25 glucose, 0.6 ascorbic acid, 3.1 pyruvic acid (bubbled with 95% O₂ and 5% CO₂, pH 7.4). Acute horizontal [for patch recordings of medial entorhinal cortex (mEC) layer II stellate neurons] or coronal slices [for patch recordings of striatal medium spiny neuron (MSNs)] at 300- μ m thickness were cut on a vibratome before transferred to artificial cerebrospinal fluid (ACSF) containing (in mmol/L): 125 NaCl, 2.5 KCl, 2.0 CaCl₂, 2.0 MgCl₂, 25 NaHCO₃, 1.25 NaH₂PO₄, 10 glucose (bubbled with 95% O₂ and 5% CO₂, pH 7.4). Brain slices were incubated at 37 °C for 20–30 min and stored at room temperature before use.

All somatic whole-cell patch-clamp recordings were performed in identified mEC layer II stellate neurons or striatal MSNs. The selection criteria for stellate cells were based on their morphological characteristics⁴⁴ (*i.e.*, large cell body presenting polygon or rhombus) and firing properties and shape of APs (*i.e.*, upon subthreshold depolarization and prominent membrane potential sags induced by both hyperpolarizing and depolarizing current injection at the somata). The selection criteria for MSNs were based on morphological characteristics with a large cell body presenting polygon or rhombus and numerous dendritic spines and their much lower resting membrane potential (RMP, about –80 mV)⁴⁵.

For whole-cell current-clamp recordings, the internal solution contained (in mmol/L): 118 KMeSO₄, 15 KCl, 2 MgCl₂, 0.2 EGTA, 10 4-(2-hydroxyethyl)-1-piperazineethanesulfonic acid (HEPES), 4 Na₂ATP, 0.3 Tris-GTP, 14 Tris-phosphocreatine, adjusted to pH 7.3 with KOH. AP5 (NMDA receptor antagonist, 50 μ mol/L), CNQX (AMPA/kainate glutamate receptor antagonist, 10 μ mol/L) were applied to block excitatory synaptic transmission, bicuculline (GABA_A receptor antagonist, 10 μ mol/L) and CGP55845 [selective γ -aminobutyric acid B (GABA_B) receptor antagonist, 2 μ mol/L] were applied to block inhibitory synaptic transmission.

The input resistance (IR) was calculated with Eq. (2):

$$\text{Input resistance} = (V_{\text{baseline}} - V_{\text{steady-state}}) \times 10 (\text{M}\Omega) \quad (2)$$

where V_{baseline} is the resting membrane potential or –70 mV (for stellate cells) or –80 mV (for MSNs) and $V_{\text{steady-state}}$ is the voltage recorded at 0–10 ms before the end of –100 pA stimulus.

For whole-cell voltage-clamp recordings of miniature inhibitory postsynaptic currents (mIPSCs), the internal solution contained (in mmol/L): 122 CsCl, 1 CaCl₂, 5 MgCl₂, 10 EGTA, 10 HEPES, 4 Na₂ATP, 0.3 Tris-GTP, 14 Tris-phosphocreatine, adjusted to pH 7.3 with CsOH. TTX (sodium channel blocker, 0.5 μ mol/L), AP5 (NMDA receptor antagonist, 50 μ mol/L), CNQX (AMPA/kainate glutamate receptor antagonist, 10 μ mol/L) for block of excitatory synaptic transmission. For

recordings of miniature excitatory postsynaptic currents (mEPSCs), the internal solution contained (in mmol/L): 118 KMeSO₄, 15 KCl, 2 MgCl₂, 0.2 EGTA, 10 HEPES, 4 Na₂ATP, 0.3 Tris-GTP, 14 Tris-phosphocreatinin, adjusted to pH 7.3 with KOH. TTX (sodium channel blocker, 0.5 μmol/L), bicuculline (GABA_A receptor antagonist, 10 μmol/L) and CGP55845 (selective GABA_B receptor antagonist, 2 μmol/L) were applied to block inhibitory synaptic transmission. Only a single neuron was recorded in each slice.

Thick-wall borosilicate pipettes were used with open-tip resistance of 4–6 MΩ. All recordings were started at least 10 min after breakthrough for internal solution exchange equilibrium using a MultiClamp 700B amplifier (Molecular Device). Data were acquired with a sampling rate at 33 kHz with a Digidata 1440A digitizer (Molecular Devices) before filtered at 2 kHz and analyzed using pCLAMP 10.6 software. Slices were maintained under continuous perfusion of artificial cerebrospinal fluid (ACSF) at 36.5–37.5 °C with a 2–3 mL/min flow. In the whole-cell configuration, the series resistance (R_s) was 15–30 MΩ, and recordings with unstable R_s or a change of R_s > 20% were aborted.

For cell labeling, the internal solution either for whole-cell current-clamp recordings or for voltage-clamp recordings contained 0.1%–0.2% (w/v) neurobiotin tracer. At the end of electrophysiological recordings (~30 min), slices were stored in 4% paraformaldehyde overnight at 4 °C before processed with streptavidin conjugated to Alexa 488 (1:500; Invitrogen) or Alexa 594 (1:500; Invitrogen) for neurobiotin visualization⁴⁶.

2.14. Imaging and Sholl analysis

Neurobiotin labeled neurons in brain slices were imaged using a laser scanning confocal microscope (Leica Microsystems) with ×63 oil-immersion objectives for stellate neurons or MSNs and ×40 oil-immersion objectives for cortical layer V pyramidal neurons. For measurement of total dendritic branch length, 3D Z-stacks were reconstructed from ~40–250 sequential scans at low zoom (1024 × 1024 μm) at 1-μm steps. The concentric sphere method of Sholl⁴⁷ was used to analyze dendritic complexity. Briefly, concentric circles at 20 μm intervals were drawn around the somata and the number of dendrites crossing each circle was counted manually from the two-dimensional display. Dendritic spine images were captured with ×63 oil-immersion objectives at 0.3-μm steps before further 6× high zoom amplification. Spine density was sampled in second- to fourth-order branches. The dendrite derived directly from the soma is the first-order branch or primary dendrite. The daughter branches arising from the first-order branch are second-order branches, and so on. Results were expressed in terms of spine density/10 μm.

2.16. Multi-electrode array (MEA) recordings of primary cortical neurons

Cortical neurons from newborn C57BL/6J mice (< 24 h) were acutely dissociated and suspended in NbActiv4 (Brain Bits) medium. Approximately $\sim 7 \times 10^4$ neurons in 8 μL medium were seeded on 12-well MEA plates (Axion Biosystems) coated with poly-D-lysine (40 $\mu\text{g}/\text{mL}$)/laminin (20 $\mu\text{g}/\text{mL}$). The cells were incubated at 37 °C in 5% CO_2 for 1 h. NbActiv4 media (0.4 mL per well) were added to the cells. On the 3rd day, cytarabine (2.5 $\mu\text{g}/\text{mL}$) was added to suppress the proliferation of glial cells by mitotic inhibition for up to seven days⁴⁸. Cells were maintained at 37 °C in a 5% CO_2 incubator, and culture medium was changed every 3 days by replacing approximately 2/3 of the medium.

Neurons were in fresh culture medium for 1 h before recorded in a 12-well plate consisting of a total of 768 electrodes. We recorded 30 min before and record 60 min after FB (100 $\mu\text{mol}/\text{L}$ in neuronal medium) application using an MEA system (Axion Biosystems). For each experiment, at least 7–13 wells (with 448–832 available electrodes for recording) were used to assess neuronal excitability. A spike detection criterion of >6 standard deviations above background signals were used to separate monophasic or biphasic action potential spikes from noise. Active electrodes were defined as 1 spike over a 200 s analysis period⁴⁹. Firing frequencies were averaged among all active electrodes from wells. MEA data were analyzed using Axion Integrated Studio AxIS2.1 (Axion Biosystems) and NeuroExplorer (Nex Technologies). The recording files were divided into every 5-min. The average response for each 5-min was calculated, and the lowest value was selected for analysis. The neuronal and network activities were assessed by criteria such as spike frequency, burst frequency, network burst frequency and synchrony index. A minimum of 5 spontaneous spikes with a maximum inter-spike interval of 0.1 s was defined as a burst, as previously reported⁵⁰, and a minimum of 10 spontaneous spikes with a maximum inter-spike interval of 0.1 s and a minimum of 25% active electrodes were as a network burst. Specifically, the network burst was manually verified by a single-blind method.

2.17. Statistical analysis

Unless otherwise noted, all data were expressed as the means \pm standard error of mean (SEM), statistical significance was determined using unpaired Student's *t*-test for the comparison between two groups. Two-way ANOVA was used for the comparison between multiple groups. Each “*n*” indicates an independent experiment, and $P < 0.05$ was statistically significant.

3. Results

3.1. Upregulation of TRPV3 channel expression in mouse model of cerebral ischemia/reperfusion injury

To recapitulate the cerebral expression of TRPV3^{25,26,51-53}, we started examining TRPV3 channel proteins in mouse cerebral neurons using immunohistochemistry. As shown in Fig. 1A, TRPV3 channel proteins (green) were robustly expressed in cultured mouse cortical neurons (red) as detected by TRPV3 specific monoclonal antibodies that were also used for detection of mTRPV3 proteins expressed in HEK 293 cells in immunofluorescence assay (Supporting Information Fig. S1A), confirming the neuronal expression of TRPV3 channel proteins.

To examine TRPV3 expression after ischemia, we established the mouse model of ischemia/reperfusion (I/R) injury induced by transient middle cerebral artery occlusion (tMCAO) under the monitoring of local cerebral blood flow (LCBF) by laser Doppler flowmetry during surgery with 80% decline of LCBF over baseline in either WT *Trpv3*^{+/+} and *Trpv3*^{-/-} (pore-deleted mutant) mice (Fig. 1B). The *Trpv3* mRNA level was increased in the ischemic regions at 6 and 12 h after 1.5 h-tMCAO injury (Fig. 1C and Fig. S1B). TRPV3 protein expressions were also increased by 1.8-fold at 6 h, 2.3-fold at 12 h and 4.9-fold at 24 h (Fig. 1D) or 7.1-fold in the penumbra and 14.5-fold in the ischemic core as compared with the contralateral regions at 24 h reperfusion (Fig. 1E). Cerebral I/R injury also resulted in an upregulation of truncated TRPV3 proteins in the ischemic core and penumbra in *Trpv3*^{-/-} brain (Fig. 1F), consistent with the observations from WT mice (Fig. 1E). These results indicate a time-dependent upregulation of TRPV3 induced by cerebral I/R injury, suggesting a role of TRPV3 channel in progressive neuronal death after ischemic stroke.

Insert Fig. 1

3.2. Attenuation of ischemic injury in *Trpv3*^{-/-} mice

To examine the functional role of TRPV3 in ischemic injury, we utilized *Trpv3* null (*Trpv3*^{-/-}) mice in which the 13 and 14 exons of *Trpv3* gene encoding the pore region were genetically deleted by expressing the targeted knockout construct²⁸. Homozygous *Trpv3*^{-/-} mice expressed a low level of truncated TRPV3 transcripts without pore region detected by TRPV3 monoclonal antibody recognizing the antigen of residues 458–474 between S1 to S2 transmembrane domain (Fig. 2A). Neurological scoring revealed that *Trpv3*^{-/-} mice exhibited a significant reduction of neurological deficit scores on Day 1, 3 and 5 after tMCAO (Fig. 2B). The brain infarct volume (8.7±2.6%) was significantly reduced in *Trpv3*^{-/-} mice as compared with *Trpv3*^{+/+} mice (31±4.9%, Fig. 2C), and the cerebral edema in *Trpv3*^{-/-} mice was also significantly attenuated with brain water content at

79.6±0.3%, as compared with 81.4±0.5% in *Trpv3*^{+/+} mice (Fig. 2D). These *in vivo* results demonstrated that *Trpv3* deficiency could protect against ischemic injury.

As prolonged oxygen–glucose deprivation (OGD) can also cause neurotoxicity, we further tested the effect of TRPV3 silencing on cortical neurons subjected to OGD insult for 30 min and reoxygenation for 24 h. As shown in Fig. 2E–G, increase of lactate dehydrogenase (LDH) release and reduction of cell viability were observed in three different types of neurons subjected to OGD/R damage. In contrast to *Trpv3*^{+/+} neurons, these detrimental effects were significantly reduced in *Trpv3*^{+/-} and *Trpv3*^{-/-} neurons, further demonstrating that suppressing TRPV3 channel function attenuated OGD-induced neurotoxicity.

Insert Fig. 2

3.3. Reduction of neuronal excitability and excitatory transmissions of striatal medium spiny neurons and cortical stellate neurons by *Trpv3* silencing

As cortical neurons surrounding acute ischemic infarcts undergo repetitive spontaneous depolarizations⁵⁴, we reasoned that silencing of TRPV3 might affect neuronal excitability. To test this notion, we carried out brain-slice recordings of principal medium spiny neurons (MSNs, Fig. 3A) from the striatum which is vulnerable to hypoxia and ischemia⁵⁵. In response to a family of 400 ms current steps, MSNs from *Trpv3*^{-/-} mice fired significantly fewer action potentials (Aps, Fig. 3B–C), and exhibited hyperpolarized resting membrane potential (RMP, Fig. 3D and Supporting Information Table S1). To exclude the influence of voltage-gated ion channels, we held neurons at the membrane potential of –80 mV and recorded APs that showed no difference between *Trpv3*^{-/-} and *Trpv3*^{+/+} neurons (Fig. 3E and F), demonstrating that *Trpv3*^{-/-} MSNs were less excitable due to their hyperpolarized RMP.

To further explore the impact of *Trpv3* null on synaptic transmissions, we recorded miniature excitatory postsynaptic currents (mEPSCs) and miniature inhibitory postsynaptic currents (mIPSCs) from MSNs of brain slices using voltage-clamp recordings. As shown in Fig. 3G–I and Supporting Information Table S2, the mEPSC frequency of *Trpv3*^{-/-} neurons was decreased with its cumulative probability moving towards a longer inter-event interval. While the mIPSC frequency was not change between *Trpv3*^{+/+} and *Trpv3*^{-/-} MSNs, which means the *Trpv3* null decreased the excitatory but not the inhibitory synaptic transmissions. Sholl analysis of dendritic trees further showed no observable differences in dendritic trees of stellate neurons in mEC layer II, pyramidal neurons in cortical layer V and MSNs in the striatum between *Trpv3*^{+/+} and *Trpv3*^{-/-} neurons (Supporting Information Fig. S2), suggesting that *Trpv3* null had little influence on neural development. The

dendrite spine density of *Trpv3*^{-/-} MSNs in striatum was decreased across classifications (2° and 3°, Fig. 3J), consistent with the reduced mEPSC frequency.

Insert Fig. 3

We also recorded layer II principal stellate neurons from the medial entorhinal cortex (mEC) that provides the main excitatory inputs to the hippocampus⁵⁶ (Supporting Information Fig. S3A). Current-clamp recordings of cortical neurons in *Trpv3*^{-/-} brain slices revealed less firing of APs with lower RMP and smaller input resistance (IR, Fig. S3B–S3E), indicating that cortical stellate *Trpv3*^{-/-} neurons were less excitable. The mEPSC frequency of *Trpv3*^{-/-} stellate neurons was also decreased with a longer inter-event interval cumulative probability (Fig. S3F–S3H). And the dendrite spine density was also reduced across classifications (2°, 3° and 4°, Fig. S3I), consistent with the results from *Trpv3*^{-/-} MSNs. Other kinetic parameters of mEPSCs and all kinetic parameters of mIPSC of stellate neurons or MSNs were not significantly different between the two groups (Table S2). These results demonstrate that *Trpv3* null striatal MSNs or cortical stellate neurons exhibited a reduction of both intrinsic neuronal excitability and excitatory transmissions.

3.4. Overexpressing TRPV3 increases neural excitability and aggravates ischemic injury

Earlier data show that silencing of TRPV3 reduced ischemic injury and neural excitability, we wondered if overexpressing TRPV3 could lead to an opposite effect. To examine the impact of TRPV3 overexpression on neural excitability, we carried out bilateral intracerebral injections of adeno-associated virus (AAV) constructs of control-EGFP (AAV-Ctrl-EGFP) or TRPV3-EGFP (AAV-TRPV3-EGFP) into mouse cortical and striatal regions (Fig. 4A). Both Ctrl-EGFP and TRPV3-EGFP proteins were robustly expressed in both cortex and striatum 30 days post-injection (Fig. 4Bi and Ci). Fluorescence imaging reveals that the control-EGFP proteins only were expressed in neuronal somata (Fig. 4Bii and 4Biii), whereas TRPV3-EGFP proteins were expressed in neuronal somata, dendrites and dendritic spines (Fig. 4Cii, 4Ciii and 4Civ). TRPV3 channel overexpression in the mouse cortex or striatum was further confirmed by Western blot and real-time PCR analysis (Supporting Information Fig. S4A–S4F). Current-clamp recordings of striatal MSNs overexpressing TRPV3 channels exhibited more firing of APs with depolarized RMP (Fig. 4D–G and Supporting Information Table S3), demonstrating that MSNs overexpressing TRPV3 were hyperexcitable.

Consistent with the observations of intrinsic hyperexcitability, overexpressing TRPV3 channels also resulted in an increased mEPSC frequency (Fig. 4H–J and Table S4), and upregulation of dendrite spine density (Fig. 4K) without significant changes of dendritic tree (Fig. 4G). In the cerebral I/R injury *in vivo* model, overexpressing TRPV3 exacerbated the neurological deficit scores,

increased the brain water content and also the infarct size (Fig. 4L–N). These results indicate that overexpressing TRPV3 caused neural hyperexcitability and aggravated ischemic injury.

Insert Fig. 4

3.5. Re-expressing TRPV3 reverses neural excitability and neuroprotection against cerebral I/R injury

To further confirm the causative role of TRPV3 in the cerebral I/R injury, we re-expressed TRPV3 channels in *Trpv3* null mice by intracerebral injections of AAV-TRPV3-EGFP or AAV-Ctrl-EGFP constructs into left cortical and striatal regions because we carried out the left MCAO model (Fig. 5A). Current-clamp recordings of MSNs from *Trpv3* gene knock-in (*Trpv3*^{-/-} + TRPV3) mice showed an increased firing, depolarized RMP and increased input resistance as compared with *Trpv3* KO neurons infected with AAV-Ctrl (*Trpv3*^{-/-} + Ctrl) (Fig. 5B–F and Supporting Information Table S5), indicating that genetic rescue by re-expressing TRPV3 increased neuronal excitability of *Trpv3* null neurons.

Re-expressing TRPV3 also resulted in an increased mEPSC frequency (Fig. 5G–I and Supporting Information Table S6) and augmented dendrite spine density (Fig. 5J) without dendritic tree change (Fig. S4H), demonstrating that re-expression of TRPV3 channels rescued the decreased excitatory network excitability and dendrite spine density of *Trpv3* null MSNs. *In vivo* evaluation further revealed that re-expressing TRPV3 also significantly exacerbated the neurological deficit, increased the brain water content and the infarct area (Fig. 5K–M), further confirming that re-expressing TRPV3 channels in *Trpv3* null mouse brain reinstated the susceptibility to ischemic damage.

Insert Fig. 5

3.6. Pharmacological inhibition of TRPV3 channels reduces neural excitability and alleviates cerebral I/R injury in mice

To examine the effect of pharmacological inhibition of TRPV3 on neural excitability, we recently identified a natural TRPV3 inhibitor forsythoside B from plant *Lamiophlomis rotate* that specifically inhibits TRPV3 current activated by 2-aminoethoxydiphenyl borate (2-APB) at 50 $\mu\text{mol/L}$ with an IC_{50} of 6.7 $\mu\text{mol/L}$ without obvious inhibitory effects on other thermoTRP channels such as hTRPA1, hTRPV1 and hTRPV4⁵⁷. We recorded the electrophysiological signals of mouse primary cortical neurons in the presence of forsythoside B using the multi-electrode array (MEA) assay. Application of forsythoside B (100 $\mu\text{mol/L}$) resulted in a decrease of neuronal firing, burst frequency and

synchrony index (Fig. 6A, B, and Supporting Information Fig. S5A–S5D). Brain slice recordings further confirmed that perfusion of forsythoside B (50 $\mu\text{mol/L}$) decreased the mEPSC frequency and the RMP in WT + Ctrl and WT + TRPV3 infected MSNs (Fig. 6C–F, Supporting Information Tables S7 and S8), meanwhile FB also decreased the mEPSC frequency in WT stellate neurons (Fig. S5E–S5H).

We also tested the effect of TRPV3 inhibition on primary cortical neurons in the presence of different concentrations of FB (0.1, 1, and 10 $\mu\text{mol/L}$) added to the culture medium 30 min before exposure to OGD for 20 min and reoxygenation of 24 h. As shown in Fig. 6G, adding forsythoside B or antioxidant edaravone (10 $\mu\text{mol/L}$) as positive control significantly increased the number of viable cells subjected to OGD injury in CCK-8 assay as compared with OGD vehicle group, demonstrating the protective effect of TRPV3 inhibition on OGD-induced neuronal injury.

We next made intravenous injections of different concentrations of forsythoside B (10, 20 and 40 mg/kg) or edaravone (3 mg/kg) at the beginning of reperfusion after mouse tMCAO for 1.5 h. Systematic administration of forsythoside B resulted in a significant reduction of infarct volumes and neurological scores after 24 h (Fig. 6H–J), but no additive effect on *Trpv3* null mice (Fig. 6K–M). All these results indicate that pharmacological inhibition of TRPV3 reduced neural hyperexcitability and attenuated ischemic brain injury.

To determine the therapeutic time window of FB, we tested the time-dependent effect of FB on the viability of primary mouse cortical neurons subjected to OGD/R in CCK8 assay. As shown in supplementary Fig. 6, addition of FB (10 $\mu\text{mol/L}$) led to a significant alleviation of neuronal OGD/R injury at time points of 0.5 and 1 h, but not after 2 and 6 h. These results suggest that inhibition of TRPV3 can alleviate neuronal cell death after an early onset of OGD injury.

Insert Fig. 6

4. Discussion

An excessive rise in intracellular Ca^{2+} during ischemia stroke has previously been shown to be attributed to the NMDA receptor mediated current and excitotoxicity⁵⁸. However, blocking glutamate receptors has failed to translate any clinical benefits for the last three decades as the high glutamate concentration has a narrow window of only one hour post-stroke⁵⁹, suggesting a sustained Ca^{2+} entry after ischemic injury is likely mediated by a non-NMDA mechanism.

In this study we attempted to identify and validate molecular targets underlying the sustained calcium entry that causes progressive brain cell death after stroke. Our rationale is based on the observations that ischemic stroke results in acute brain tissue acidosis due to elevated lactic acid in

the affected area. Lactic acid, an alpha hydroxyl acid, is constantly produced from pyruvate by lactate dehydrogenase during normal metabolism; however, its production dramatically increases as ischemia suppresses oxygen-dependent acetyl CoA synthesis process. Although the acetyl CoA-based citric acid cycle is usually the main energy source for living tissues, the latest findings indicate that in mammalian brain lactate is preferentially metabolized⁶⁰⁻⁶³. The resultant strenuous anaerobic process causes acute brain acidosis⁶⁴. Lactic acid-induced acidosis occurring in well-oxygenated muscles during intense exercise can increase blood lactic acid concentration from 1–2 to over 20 mmol/L. In blood supply-deprived brain regions, lactic acid synthesis elevates while its removal is absent. Therefore, lactic acid accumulation can be even more severe. Indeed, upon ischemia onset, within minutes both extracellular and intracellular pH of the affected region can drop below 6.5, and accumulation of CO₂ following ischemic stroke also contributes to a local decrease in pH^{65,66}.

The warm-temperature activated Ca²⁺-permeable TRPV3 channel expressed in the brain can be directly activated by intracellular acidification^{24,67}, suggesting that TRPV3 can be critical to induce cytotoxicity and cerebral tissue damage after ischemic stroke (Fig. 7). Genetic gain-of-functional mutations of TRPV3 lead to apoptosis of keratinocytes and skin diseases in patients with Olmsted syndrome⁶⁸, although there is no clinical evidence yet for their susceptibility for ischemic stroke. In hippocampal neurons, TRPV3 has been indicated to contribute to ischemia-induced depolarization, intracellular Ca²⁺ accumulation and cell swelling⁵². When hippocampal slices are cooled slightly or when ruthenium red was applied to block TRPV3, these harmful cellular effects are sharply reduced⁵². Indeed, before TRPV3 and other heat-sensitive TRP channels were discovered, a number of independent studies show that mild hypothermia can protect CA1 neurons from damage caused by experimental ischemia⁶⁹⁻⁷⁵. These observations may help explain the beneficial effects of deliberate hypothermia treatment of patients with head injuries^{72,76}. ThermoTRPV3 is broadly distributed in brain tissues⁷⁷, and also is present throughout the cortex, thalamus, striatum and cerebellum^{25,26,51,52}, among which its expression and function in hippocampal pyramidal neurons are studied^{25,26,51,78}. Pyramidal neurons lacking thermoTRPV1 or TRPV4 are less excitable, with resting membrane potentials about 5 mV more negative than cells from WT mice^{79,80}. Application of ruthenium red, a TRPV channel blocker, hyperpolarizes the membrane of WT cells, further supporting that these TRPV channels are open at the resting membrane potential and contributing to depolarization of brain neurons⁵².

TRPV3 channel is activated by arachidonic acid metabolites such as epoxyeicosatrienoic acids that present in cerebral neurons^{81,82}, and can be potentiated by free fatty acids⁸³. The basal activity of

TRPV3 makes neurons more susceptible to acidification attack because the channel can conduct H^+ current to cause intracellular acidification. The high susceptibility of TRPV3 to a variety of physical and chemical stimuli makes it a very attractive target for clinical intervention under ischemia conditions. It has been reported that a natural incensole acetate, a *Boswellia* resin constituent, shows protective effects on cerebral ischemic injury⁸⁴. The incensole acetate, as a non-specific TRPV3 agonist, activates TRPV3 expressed in HEK293 cells at high concentration of 200 $\mu\text{mol/L}$ ⁸⁴, and also inhibits inflammatory mediators, including $\text{TNF-}\alpha$, $\text{IL-1}\beta$, $\text{TGF-}\beta$ and $\text{NF-}\kappa\text{B}$ to reduce inflammation in cerebral ischemic injury⁸⁵.

While most H^+ -activated receptors and channels in the brain can sense extracellular acidification, the warm-temperature sensitive and Ca^{2+} -permeable TRPV3 is an only known member of TRPs activated by intracellular protons²⁴. This distinct feature makes neuronal TRPV3 an interesting and attractive target for stroke pathology and recovery as well as potential intervention. In support of this view, a low extracellular pH under aerobic conditions has little impact on intracellular calcium $[\text{Ca}^{2+}]_i$, whereas hypoxia significantly increases Ca^{2+} influx in neurons and decreases extracellular pH⁸⁶. These observations define TRPV3 as a unique brain pH-sensitive channel that mediates Ca^{2+} influx. As a non-selective cation channel, TRPV3 is featured of high PCa/PNa ratio of 3-to-12^{25,27}. Importantly, TRPV3 does not desensitize, and instead, repetitive or prolonged stimulation causes strong potentiation of TRPV3 activation, making the Ca^{2+} -carrying current larger over time^{87,88}. In contrast, most ligand-gated ion channels, including NMDA receptors, can rapidly desensitize upon the persistent presence of ligand molecules, often at a time constant in the range of seconds to minutes.

5. Conclusions

We propose that overactive intracellular proton-sensitive and calcium-permeable TRPV3 plays a causative role in neuronal hyperexcitability and ischemia-induced sustained Ca^{2+} influx for cerebral injury after stroke (Fig. 7), a mechanism that is independent of NMDA-mediated excitotoxicity. Unlike most other neurological behavioral tests, the neurological deficit scores were determined by Longa's method³², which is positively related to neuronal death⁸⁹, rather than directly to the neuronal excitabilities. Our data indicate that the over-expression of TRPV3 channels increases neuronal excitabilities, resulting in those neurons more vulnerable to excitatory toxicity or death due to ischemic injury and leading to more severe neurological deficit scores. In consideration of thermoTRPV3 as an attractive drug target with therapeutic potential, we further screened out and identified a natural compound forsythoside B that reduces neuronal hyperexcitability and alleviates

ischemic injury. Therefore, pharmacological inhibition of overactive TRPV3 after ischemic injury may lead to an effective prevention and therapy for ischemic brain stroke.

Acknowledgments

This work was supported by grants awarded to K.W. from National Natural Science Foundation of China (81903734 and 81973299) and the Ministry of Science and Technology of China (2018ZX09711001-004-006).

Author contributions

Xiaoling Chen designed and performed experiments of immunostaining, Western blot, MCAO model, OGD model, MEA recording, dendritic spine, and drafted the manuscript. Jingliang Zhang designed and performed brain slice electrophysiological recordings, and drafted the manuscript. KeWei Wang designed the study and finalized the manuscript. All authors approved the final version of the manuscript.

Conflicts of interest

The authors declare no competing interests.

References

- 1 Rehncrona S. Brain acidosis. *Ann Emerg Med* 1985; **14**: 770-6.
- 2 Simon R, Xiong Z. Acidotoxicity in brain ischaemia. *Biochem Soc Trans* 2006; **34**: 1356-61.
- 3 Huang Y, McNamara JO. Ischemic stroke: "acidotoxicity" is a perpetrator. *Cell* 2004; **118**: 665-6.
- 4 Chu XP, Xiong ZG. Acid-sensing ion channels in pathological conditions. *Adv Exp Med Biol* 2013; **961**: 419-31.
- 5 Siesjo BK, Katsura K, Kristian T. Acidosis-related damage. *Adv Neurol* 1996; **71**: 209-36.
- 6 Wang WZ, Chu XP, Li MH, Seeds J, Simon RP, Xiong ZG. Modulation of acid-sensing ion channel currents, acid-induced increase of intracellular Ca^{2+} , and acidosis-mediated neuronal injury by intracellular pH. *J Biol Chem* 2006; **281**: 29369-78.
- 7 Guo Y, Zhou IY, Chan ST, Wang Y, Mandeville ET, Igarashi T, et al. pH-sensitive MRI demarcates graded tissue acidification during acute stroke—pH specificity enhancement with magnetization transfer and relaxation-normalized amide proton transfer (APT) MRI.

- Neuroimage* 2016; **141**: 242-9.
- 8 Xiong ZG, Zhu XM, Chu XP, Minami M, Hey J, Wei WL, et al. Neuroprotection in ischemia: blocking calcium-permeable acid-sensing ion channels. *Cell* 2004; **118**: 687-98.
- 9 Xiong ZG, Chu XP, Simon RP. Acid sensing ion channels—novel therapeutic targets for ischemic brain injury. *Front Biosci* 2007; **12**: 1376-86.
- 10 Xiong ZG, Pignataro G, Li M, Chang SY, Simon RP. Acid-sensing ion channels (ASICs) as pharmacological targets for neurodegenerative diseases. *Curr Opin Pharmacol* 2008; **8**: 25-32.
- 11 Nakamura K, Kamouchi M, Arimura K, Nishimura A, Kuroda J, Ishitsuka K, et al. Extracellular acidification activates cAMP responsive element binding protein via Na^+/H^+ exchanger isoform 1-mediated Ca^{2+} oscillation in central nervous system pericytes. *Arterioscler Thromb Vasc Biol* 2012; **32**: 2670-7.
- 12 Begum G, Song S, Wang S, Zhao H, Bhuiyan MIH, Li E, et al. Selective knockout of astrocytic Na^+/H^+ exchanger isoform 1 reduces astrogliosis, BBB damage, infarction, and improves neurological function after ischemic stroke. *Glia* 2018; **66**: 126-44.
- 13 Choi DW. Glutamate neurotoxicity and diseases of the nervous-system. *Neuron* 1988; **1**: 623-34.
- 14 Choi DW. Calcium-mediated neurotoxicity—relationship to specific channel types and role in ischemic damage. *Trends Neurosci* 1988; **11**: 465-69.
- 15 Lee JM, Zipfel GJ, Choi DW. The changing landscape of ischaemic brain injury mechanisms. *Nature* 1999; **399**: A7-14.
- 16 Chamorro A, Dirnagl U, Urra X, Planas AM. Neuroprotection in acute stroke: targeting excitotoxicity, oxidative and nitrosative stress, and inflammation. *Lancet Neurol* 2016; **15**: 869-81.
- 17 Ikonomidou C, Turski L. Why did NMDA receptor antagonists fail clinical trials for stroke and traumatic brain injury?. *Lancet Neurology* 2002; **1**: 383-86.
- 18 Albeni BC, Igoechi C, Janigro D, Ilkanich E. Why do many NMDA antagonists fail, while others are safe and effective at blocking excitotoxicity associated with dementia and acute injury?. *Am J Alzheimers Dis Other Demen* 2004; **19**: 269-74.
- 19 MacDonald JF, Jackson MF. Transient receptor potential channels of the melastatin family and ischemic responses of central neurons. *Stroke* 2007; **38**: 665-9.
- 20 Lai TW, Zhang S, Wang YT. Excitotoxicity and stroke: identifying novel targets for neuroprotection. *Prog Neurobiol* 2014; **115**: 157-88.

- 21 Simard JM, Tarasov KV, Gerzanich V. Non-selective cation channels, transient receptor potential channels and ischemic stroke. *Biochim Biophys Acta* 2007; **1772**: 947-57.
- 22 Zhang E, Liao P. Brain transient receptor potential channels and stroke. *J Neurosci Res* 2015; **93**: 1165-83.
- 23 Leng T, Shi Y, Xiong ZG, Sun D. Proton-sensitive cation channels and ion exchangers in ischemic brain injury: new therapeutic targets for stroke?. *Prog Neurobiol* 2014; **115**: 189-209.
- 24 Cao X, Yang F, Zheng J, Wang K. Intracellular proton-mediated activation of TRPV3 channels accounts for the exfoliation effect of alpha-hydroxyl acids on keratinocytes. *J Biol Chem* 2012; **287**: 25905-16.
- 25 Xu H, Ramsey IS, Kotecha SA, Moran MM, Chong JA, Lawson D, et al. TRPV3 is a calcium-permeable temperature-sensitive cation channel. *Nature* 2002; **418**: 181-6.
- 26 Smith GD, Gunthorpe MJ, Kelsell RE, Hayes PD, Reilly P, Facer P, et al. TRPV3 is a temperature-sensitive vanilloid receptor-like protein. *Nature* 2002; **418**: 186-90.
- 27 Peier AM, Reeve AJ, Andersson DA, Moqrich A, Earley TJ, Hergarden AC, et al. A heat-sensitive TRP channel expressed in keratinocytes. *Science* 2002; **296**: 2046-9.
- 28 Moqrich A, Hwang SW, Earley TJ, Petrus MJ, Murray AN, Spencer KS, et al. Impaired thermosensation in mice lacking TRPV3, a heat and camphor sensor in the skin. *Science* 2005; **307**: 1468-72.
- 29 Schmitteckert EM, Prokop CM, Hedrich HJ. DNA detection in hair of transgenic mice—a simple technique minimizing the distress on the animals. *Lab Anim* 1999; **33**: 385-9.
- 30 Rousselet E, Kriz J, Seidah NG. Mouse model of intraluminal MCAO: cerebral infarct evaluation by cresyl violet staining. *J Vis Exp* 2012;(69):4038.
- 31 Liu F, McCullough LD. The middle cerebral artery occlusion model of transient focal cerebral ischemia. *Methods Mol Biol* 2014; **1135**: 81-93.
- 32 Longa EZ, Weinstein PR, Carlson S, Cummins R. Reversible middle cerebral artery occlusion without craniectomy in rats. *Stroke* 1989; **20**: 84-91.
- 33 Bu Q, Liu X, Zhu Y, Liu Y, Wang Y. w007B protects brain against ischemia–reperfusion injury in rats through inhibiting inflammation, apoptosis and autophagy. *Brain Res* 2014; **1558**: 100-08.
- 34 Chiang T, Messing RO, Chou WH. Mouse model of middle cerebral artery occlusion. *J Vis Exp* 2011;(48):2761.
- 35 Jackman K, Kunz A, Iadecola C. Modeling focal cerebral ischemia *in vivo*. *Methods Mol Biol*

- 2011;**793**:195-209.
- 36 Yan H, Zhang X, Hu W, Ma J, Hou W, Zhang X, et al. Histamine H3 receptors aggravate cerebral ischaemic injury by histamine-independent mechanisms. *Nat Commun* 2014; **5**: 3334.
- 37 Chao X, Zhou J, Chen T, Liu W, Dong W, Qu Y, et al. Neuroprotective effect of osthole against acute ischemic stroke on middle cerebral ischemia occlusion in rats. *Brain Res* 2010; **1363**: 206-11.
- 38 Nunez J. Primary culture of hippocampal neurons from P0 newborn rats. *J Vis Exp* 2008;(19):895.
- 39 Hu Z, Bian X, Liu X, Zhu Y, Zhang X, Chen S, et al. Honokiol protects brain against ischemia–reperfusion injury in rats through disrupting PSD95–nNOS interaction. *Brain Res* 2013; **1491**: 204-12.
- 40 Guan L, Song Y, Gao J, Gao J, Wang K. Inhibition of calcium-activated chloride channel ANO1 suppresses proliferation and induces apoptosis of epithelium originated cancer cells. *Oncotarget* 2016; **7**: 78619-30.
- 41 Livak KJ, Schmittgen TD. Analysis of relative gene expression data using real-time quantitative PCR and the $2^{-\Delta\Delta CT}$ Method. *Methods* 2001; **25**: 402-8.
- 42 Ashwal S, Tone B, Tian HR, Cole DJ, Pearce WJ. Core and penumbral nitric oxide synthase activity during cerebral ischemia and reperfusion. *Stroke* 1998; **29**: 1037-46; discussion 47.
- 43 Memezawa H, Minamisawa H, Smith ML, Siesjo BK. Ischemic penumbra in a model of reversible middle cerebral artery occlusion in the rat. *Exp Brain Res* 1992; **89**: 67-78.
- 44 Alessi C, Raspanti A, Magistretti J. Two distinct types of depolarizing afterpotentials are differentially expressed in stellate and pyramidal-like neurons of entorhinal-cortex layer II. *Hippocampus* 2016; **26**: 380-404.
- 45 Torres-Garcia ME, Solis O, Patricio A, Rodriguez-Moreno A, Camacho-Abrego I, Limon ID, et al. Dendritic morphology changes in neurons from the prefrontal cortex, hippocampus and nucleus accumbens in rats after lesion of the thalamic reticular nucleus. *Neuroscience* 2012; **223**: 429-38.
- 46 Zhang J, Chen X, Kårbø M, Zhao Y, An L, Wang R, et al. Anticonvulsant effect of dipropofol by enhancing native GABA currents in cortical neurons in mice. *J Neurophysiol* 2018; **120**: 1404-14.
- 47 Sholl DA. Dendritic organization in the neurons of the visual and motor cortices of the cat. *J Anat* 1953; **87**: 387-406.

- 48 Schwieger J, Esser KH, Lenarz T, Scheper V. Establishment of a long-term spiral ganglion neuron culture with reduced glial cell number: effects of AraC on cell composition and neurons. *J Neurosci Methods* 2016; **268**: 106-16.
- 49 Yang Y, Adi T, Effraim PR, Chen L, Dib-Hajj SD, Waxman SG. Reverse pharmacogenomics: carbamazepine normalizes activation and attenuates thermal hyperexcitability of sensory neurons due to Nav 1.7 mutation I234T. *Br J Pharmacol* 2018; **175**: 2261-71.
- 50 Liu DC, Seimetz J, Lee KY, Kalsotra A, Chung HJ, Lu H, et al. Mdm2 mediates FMRP- and Gp1 mGluR-dependent protein translation and neural network activity. *Hum Mol Genet* 2017; **26**: 3895-908.
- 51 Kunert-Keil C, Bisping F, Krüger J, Brinkmeier H. Tissue-specific expression of TRP channel genes in the mouse and its variation in three different mouse strains. *BMC genomics* 2006; **7**: 1-14.
- 52 Lipski J, Park TI, Li D, Lee SC, Trevarton AJ, Chung KK, et al. Involvement of TRP-like channels in the acute ischemic response of hippocampal CA1 neurons in brain slices. *Brain Res* 2006; **1077**: 187-99.
- 53 Caterina MJ. Transient receptor potential ion channels as participants in thermosensation and thermoregulation. *Am J Physiol Regul Integr Comp Physiol* 2007; **292**: R64-76.
- 54 Nedergaard M, Hansen AJ. Characterization of cortical depolarizations evoked in focal cerebral ischemia. *J Cereb Blood Flow Metab* 1993; **13**: 568-74.
- 55 Calabresi P, Pisani A, Mercuri NB, Bernardi G. On the mechanisms underlying hypoxia-induced membrane depolarization in striatal neurons. *Brain* 1995; **118**: 1027-38.
- 56 Steward O. Topographic organization of the projections from the entorhinal area to the hippocampal formation of the rat. *J Comp Neurol* 1976; **167**: 285-314.
- 57 Zhang H, Sun X, Qi H, Ma Q, Zhou Q, Wang W, et al. Pharmacological inhibition of the temperature-sensitive and Ca²⁺-permeable transient receptor potential vanilloid TRPV3 channel by natural forsythoside B attenuates pruritus and cytotoxicity of keratinocytes. *J Pharmacol Exp Ther* 2019; **368**: 21-31.
- 58 Tu W, Xu X, Peng L, Zhong X, Zhang W, Soundarapandian MM, et al. DAPK1 interaction with NMDA receptor NR2B subunits mediates brain damage in stroke. *Cell* 2010; **140**: 222-34.
- 59 Ikonomidou C, Turski L. Why did NMDA receptor antagonists fail clinical trials for stroke and traumatic brain injury?. *Lancet Neurol* 2002; **1**: 383-6.
- 60 Zilberter Y, Zilberter T, Bregestovski P. Neuronal activity *in vitro* and the *in vivo* reality: the

- role of energy homeostasis. *Trends Pharmacol Sci* 2010; **31**: 394-401.
- 61 Gladden LB. Lactate metabolism: a new paradigm for the third millennium. *J Physiol* 2004; **558**: 5-30.
- 62 Pellerin L, Bouzier-Sore AK, Aubert A, Serres S, Merle M, Costalat R, et al. Activity-dependent regulation of energy metabolism by astrocytes: an update. *Glia* 2007; **55**: 1251-62.
- 63 Wyss MT, Jolivet R, Buck A, Magistretti PJ, Weber B. *In vivo* evidence for lactate as a neuronal energy source. *J Neurosci* 2011; **31**: 7477-85.
- 64 Siesjö B, Ekholm A, Katsura K, Theander S. Acid-base changes during complete brain ischemia. *Stroke* 1990; **21**: III194-9.
- 65 Lam TI, Wise PM, O'Donnell ME. Cerebral microvascular endothelial cell Na/H exchange: evidence for the presence of NHE1 and NHE2 isoforms and regulation by arginine vasopressin. *Am J Physiol Cell Physiol* 2009; **297**: C278-89.
- 66 Rehncrona S, Rosén I, Smith ML. Effect of different degrees of brain ischemia and tissue lactic acidosis on the short-term recovery of neurophysiologic and metabolic variables. *Exp Neurol* 1985; **87**: 458-73.
- 67 Gao L, Yang P, Qin P, Lu Y, Li X, Tian Q, et al. Selective potentiation of 2-APB-induced activation of TRPV1-3 channels by acid. *Sci Rep* 2016; **6**: 20791.
- 68 Lin Z, Chen Q, Lee M, Cao X, Zhang J, Ma D, et al. Exome sequencing reveals mutations in TRPV3 as a cause of Olmsted syndrome. *Am J Hum Genet* 2012; **90**: 558-64.
- 69 Onitsuka M, Mihara S, Inokuchi H, Shigemori M, Higashi H. Mild hypothermia protects rat hippocampal CA1 neurons from irreversible membrane dysfunction induced by experimental ischemia. *Neurosci Res* 1998; **30**: 1-6.
- 70 Busto R, Dietrich WD, Globus MYT, Ginsberg MD. Postischemic moderate hypothermia inhibits CA1 hippocampal ischemic neuronal injury. *Neurosci Lett* 1989; **101**: 299-304.
- 71 Busto R, Dietrich WD, Globus MY-T, Valdés I, Scheinberg P, Ginsberg MD. Small differences in intrischemic brain temperature critically determine the extent of ischemic neuronal injury. *J Cereb Blood Flow Metab* 1987; **7**: 729-38.
- 72 Clifton GL, Allen S, Barrodale P, Plenger P, Berry J, Koch S, et al. A phase II study of moderate hypothermia in severe brain injury. *J Neurotrauma* 1993; **10**: 263-71.
- 73 Clifton GL, Jiang JY, Lyeth BG, Jenkins LW, Hamm RJ, Hayes RL. Marked protection by moderate hypothermia after experimental traumatic brain injury. *J Cereb Blood Flow Metab* 1991; **11**: 114-21.

- 74 Resnick DK, Marion DW, Darby JM. The effect of hypothermia on the incidence of delayed traumatic intracerebral hemorrhage. *Neurosurgery* 1994; **34**: 252-56.
- 75 Bramlett HM, Green EJ, Dietrich WD, Busto R, GLOBUS MYT, Ginsberg MD. Posttraumatic brain hypothermia provides protection from sensorimotor and cognitive behavioral deficits. *J Neurotrauma* 1995; **12**:289-98.
- 76 Mellergård P. Changes in human intracerebral temperature in response to different methods of brain cooling. *Neurosurgery* 1992; **31**: 671-77.
- 77 Caterina MJ. Transient receptor potential ion channels as participants in thermosensation and thermoregulation. *Am J Physiol Regul Integr Comp Physiol* 2007; **292**:R64-76.
- 78 Mutai H, Heller S. Vertebrate and invertebrate TRPV-like mechanoreceptors. *Cell Calcium* 2003; **33**: 471-78.
- 79 Shibasaki K, Suzuki M, Mizuno A, Tominaga M. Effects of body temperature on neural activity in the hippocampus: regulation of resting membrane potentials by transient receptor potential vanilloid 4. *J Neurosci* 2007; **27**: 1566-75.
- 80 Edwards JG, Gibson HE, Jensen T, Nugent F, Walther C, Blickenstaff J, et al. A novel non- CB1/TRPV1 endocannabinoid- mediated mechanism depresses excitatory synapses on hippocampal CA1 interneurons. *Hippocampus* 2012; **22**: 209-21.
- 81 Xu H, Delling M, Jun JC, Clapham DE. Oregon, thyme and clove-derived flavors and skin sensitizers activate specific TRP channels. *Nat Neurosci* 2006; **9**: 628-35.
- 82 Fernandes J, Lorenzo IM, Andrade YN, Garcia-Elias A, Serra SA, Fernandez-Fernandez JM, et al. IP3 sensitizes TRPV4 channel to the mechano- and osmotransducing messenger 5'-6'-epoxyeicosatrienoic acid. *J Gen Physiol* 2008; **131**: i2.
- 83 Hu HZ, Xiao R, Wang C, Gao N, Colton CK, Wood JD, et al. Potentiation of TRPV3 channel function by unsaturated fatty acids. *J Cellular Physiology* 2006; **208**: 201-12.
- 84 Moussaieff A, Rimmerman N, Bregman T, Straiker A, Felder CC, Shoham S, et al. Incensole acetate, an incense component, elicits psychoactivity by activating TRPV3 channels in the brain. *FASEB J* 2008; **22**: 3024-34.
- 85 Moussaieff A, Yu J, Zhu H, Gattioni-Celli S, Shohami E, Kindy MS. Protective effects of incensole acetate on cerebral ischemic injury. *Brain Res* 2012; **1443**: 89-97.
- 86 O'Donnell BR, Bickler PE. Influence of pH on calcium influx during hypoxia in rat cortical brain slices. *Stroke* 1994; **25**: 171-77.
- 87 Chung MK, Guler AD, Caterina MJ. Biphasic currents evoked by chemical or thermal activation of the heat-gated ion channel, TRPV3. *J Biol Chem* 2005; **280**: 15928-41.

- 88 Cheng W, Yang F, Liu S, Colton CK, Wang C, Cui Y, et al. Heteromeric heat-sensitive transient receptor potential channels exhibit distinct temperature and chemical response. *J Biol Chem* 2012; **287**: 7279-88.
- 89 Senda DM, Franzin S, Mori MA, de Oliveira RM, Milani H. Acute, post-ischemic sensorimotor deficits correlate positively with infarct size but fail to predict its occurrence and magnitude after middle cerebral artery occlusion in rats. *Behav Brain Res* 2011; **216**: 29-35.
- 90 Siesjo BK, Katsura K, Mellergard P, Ekholm A, Lundgren J, Smith ML. Acidosis-related brain damage. *Progress Brain Res* 1993; **96**: 23-48.
- 91 Broad LM, Mogg AJ, Eberle E, Tolley M, Li DL, Knopp KL. TRPV3 in drug development. *Pharmaceuticals (Basel)* 2016; **9**: 55.
- 92 Nedergaard M, Kraig RP, Tanabe J, Pulsinelli WA. Dynamics of interstitial and intracellular pH in evolving brain infarct. *Am J Physiol* 1991; **260**: R581-R88.

Figures captions

Figure 1 Upregulation of TRPV3 channel expressions induced by cerebral ischemic–reperfusion injury in mice. (A) Upper panels, immunohistochemical co-staining of TRPV3 protein (green) and NeuN biomarker (red) in cultured mouse primary cortical neurons were observed using confocal microscopy with 4 repeats. Bottom panels, secondary antibodies were used as a negative control. Nuclei are stained by DAPI (blue). (B) Measurement of ipsilateral local cerebral blood flow (LCBF) by laser Doppler flowmetry in *Trpv3*^{+/+} mouse (left panel) and *Trpv3*^{-/-} mouse (middle panel) subjected to 1.5 h ischemia/ 24 h reperfusion (I/R) injury induced by transient middle cerebral artery occlusion (tMCAO) surgery. Up- and down-arrows (red) indicate the insertion and withdrawal of the filament, respectively. In right panel, the statistics of declined LCBF of baseline. Ns, no significance. (C) Top panel, a timeline for generation of tMCAO in mice subjected to 1.5 h ischemia (I) before 6, 12 or 24 h reperfusion (R) injury. Bottom left panel, representative image of ischemic (Isc) and contralateral (Con) regions at 24 h reperfusion, and upregulation of *Trpv3* mRNA levels in ischemia region at 6 and 12 h reperfusion after 1.5 h ischemia using quantitative real-time PCR assay (bottom right panel) using the two-way ANOVA followed by Holm-Sidak's test from GraphPad 8. (D) Time-dependent increase of TRPV3 protein expressions in the ischemia area at 6, 12, and 24 h of reperfusion after 1.5 h ischemia (top panel) and their analysis for TRPV3 protein expression (bottom panel) using the two-way ANOVA followed by Uncorrected Fisher's LSD from GraphPad 8. (E) Top panel, Western blot analysis for upregulation of TRPV3 protein expressions in ischemia penumbra and core regions in *Trpv3*^{+/+} mice. Bottom left panel, a representative image of ischemic core, penumbra (Pen) and contralateral (Con) regions at 24 h of reperfusion and their analysis for TRPV3 expression (bottom panel) using the two-way ANOVA followed by Uncorrected Fisher's LSD from GraphPad 8. (F) Upregulation of TRPV3 protein expressions in ischemia penumbra and core regions in *Trpv3*^{-/-} mice subjected to tMCAO injury and their analysis of TRPV3 expression using the two-way ANOVA followed by Uncorrected Fisher's LSD from GraphPad 8. Data are expressed as mean \pm SEM, *n* represents the number of mice. **P* < 0.05, ***P* < 0.01, ****P* < 0.001. Also see Supporting Information Fig. S1A for antibody specificity; Fig. S1B for upregulations of *Trpv3* mRNA levels after tMCAO injury.

Figure 2 Silencing of TRPV3 attenuates ischemic brain injury. (A) Top panel, a schematic TRPV3 topology illustrating the deletion of exons 13 and 14 that encoding the putative pore region and adjacent transmembrane segments five and six. And the region of residues 458–474 between S1 to S2 transmembrane domain recognized by the monoclonal TRPV3 antibody (ab94582). Bottom panel, the PCR genotyping for confirmation of *Trpv3*^{-/-} with detected cDNA band at 182 bp, *Trpv3*^{+/+} with 304 bp, and *Trpv3*^{+/-} mice carrying the two bands. (B) Attenuation of neurological deficit scores at time points of day 1, 3 and 5 after 1.5 h tMCAO injury in *Trpv3*^{-/-} mice. Two-tailed nonparametric Mann–Whitney test. Data are shown as the median with 95% CI, and *n* indicates the number of mice. **P* < 0.05. (C) Representative TTC-stained brain slices at 24 h after reperfusion from *Trpv3*^{+/+} and *Trpv3*^{-/-} mice. The white regions indicate the infarct size, and the red regions indicate the viable tissues. Reduction of infarction volume in *Trpv3*^{-/-} mice subjected to cerebral I/R (1.5 h/24 h) injury. (D) A decrease of brain water content in *Trpv3*^{-/-} mice subjected to cerebral I/R (1.5 h/24 h) injury. (E) Top panel, a schematic drawing for primary mouse cortical neurons subjected to 30 min oxygen-glucose deprivation and 24 h reoxygenation (OGD/R) injury. The representative images for morphological changes of *Trpv3*^{+/+} and *Trpv3*^{-/-} neurons subjected to OGD/R injury (bottom panel). Scale bar, 100 μm. The statistics of the lactate dehydrogenase (LDH) release (F) in LDH assay and cell viability (G) by cell counting kit-8 (CCK8) assay of primary cortical neurons subjected to OGD/R injury from *Trpv3*^{+/+}, *Trpv3*^{+/-} and *Trpv3*^{-/-} mice. Data are expressed as means ± SEM, *n* indicates the number of mice. Two-way ANOVA was used and followed by Dunnett's test from GraphPad 8. **P* < 0.05, ***P* < 0.01 and ****P* < 0.001.

Figure 3 Reduction of neural excitability in *Trpv3*^{-/-} striatal medium spiny neurons. (A) Morphology of a typical mouse medium spiny neuron (MSN) in the striatum labeled with neurobiotin. (B) Representative traces for neuronal firings by whole-cell current-clamp recordings of the MSN at resting membrane potential (RMP) from *Trpv3*^{+/+} and *Trpv3*^{-/-} mice. (C) The summary of action potential numbers (AP No.) at RMP. AP No. of each current step was analyzed by two-tailed nonparametric Mann–Whitney test. (D) The summary of RMP values. (E) Representative traces for neuronal firings of MSNs at -80 mV from *Trpv3*^{+/+} and *Trpv3*^{-/-} mice. (F) The summary of AP No. at -80 mV. AP No. of each current step was analyzed by two-tailed nonparametric Mann–Whitney test. (G) Representative traces for mEPSCs in striatal MSNs from *Trpv3*^{+/+} and *Trpv3*^{-/-} mice

by whole-cell voltage clamp recordings. (H) The frequency of mEPSCs and associated cumulative probability (I). (J) Left panel, representative images showing the morphology of the spines in striatal MSNs from *Trpv3^{+/+}* and *Trpv3^{-/-}* mice. The summary of dendrite spine density (middle panel) and a schematic drawing of dendrite branch classifications (right panel). Data are expressed as mean \pm SEM, *n* represents the number of neurons patched. Two-way ANOVA was used and followed by Holm–Sidak's test from GraphPad 8. **P* < 0.05, ***P* < 0.01 and ****P* < 0.001. Also see Supporting Information Fig. S2 for no obvious morphological changes of complex dendritic trees in *Trpv3^{+/+}* and *Trpv3^{-/-}* neurons; Fig. S3 for reduced neural excitability of cortical stellate neurons from *Trpv3^{-/-}* mice and Table S1 for electrophysiological parameters on medium spiny neurons (MSNs) and layer II principal stellate neurons from medial entorhinal cortex (mEC) from *Trpv3^{+/+}* and *Trpv3^{-/-}* mice and Table S2 for mEPSCs and mIPSCs parameters of mouse striatal MSNs or mEC layer II stellate neurons.

Figure 4 Overexpressing TRPV3 increases neural excitability and aggravates cerebral I/R injury. (A) Experimental timeline for viral injection on Day 0 and detection expression of AAV constructs at day 30 and executing tMCAO model on Day 31, and illustration of bilateral viral injecting sites in mouse cortex and striatum. Schematic drawing of AAV constructs for a control (AAV-Ctrl-EGFP) and overexpression of TRPV3-EGFP (AAV-TRPV3-EGFP). ITR, inverted terminal repeats; hsyn, human synapsin promoter; EGFP, enhanced green fluorescent protein. (B) Representative fluorescent images of coronal mouse brain slice are from WT mice injected with control virus after 30 days (Bi, WT + Ctrl) and control-EGFP proteins expressed in neuronal somata (Bii and Biii). The amplifications are from the sensory cortical regions indicated by white square frame. (C) Representative fluorescent coronal brain slice is from WT mice injected with TRPV3 overexpressing virus. (Ci, WT + TRPV3) and TRPV3-EGFP proteins expressed in neuronal somata (Cii), dendrites (Ciii) and dendritic spines (Civ) from Ciii. (D) Morphology of a typical mouse MSN in striatum labeled with neurobiotin that we patched. (E) Representative traces for neuronal firings in striatal MSNs from WT + Ctrl and WT + TRPV3 groups obtained at RMP by whole-cell current-clamp recordings. (F) Graph depicting the average AP No. at RMP. AP No. of each current step was analyzed by two-tailed nonparametric Mann–Whitney test. (G) The summary of RMP values. (H) Representative traces for mEPSC in striatal MSNs

from WT + Ctrl and WT + TRPV3 groups by whole-cell voltage-clamp recordings. (I) The frequency of mEPSC and associated cumulative probability (J). (K) Representative images showing the morphology of spines in striatal MSNs from WT + Ctrl and WT + TRPV3 mice. The summary of spine density and a schematic drawing of a dendrite. Two-way ANOVA followed by Holm-Sidak's test. (L) Increase of neurological deficit scores, (M) brain water content (%) and (N) infarction volume (%) in TRPV3 overexpressing mice subjected to cerebral I/R (1.5/24 h) injury. Data for neurological deficit scores are shown as the median with 95% CI, and was analyzed two-tailed nonparametric Mann–Whitney test. Data are represented as mean \pm SEM, n represents the number of neurons or the number of mice. * $P < 0.05$, ** $P < 0.01$ and *** $P < 0.001$. Also see Supporting Information Fig. S4 for efficiency of TRPV3 overexpression in mouse cortex and striatum; and Table S3 for electrophysiological parameters of mouse striatal MSNs between WT + Ctrl and WT + TRPV3 groups; Table S4 for mEPSC parameters in mouse striatal MSNs between WT + Ctrl and WT + TRPV3 groups.

Figure 5 Re-expressing TRPV3 reverses neural excitability and neuroprotection subject to cerebral I/R injury. (A) Experimental timelines for viral injections of AAV-Ctrl-EGFP and AAV-TRPV3-EGFP constructs into the left cortex and striatum in *Trpv3* null mice on Day 0 (D0), and generation of tMCAO model on Day 31 (D31). (B) Morphology of a typical mouse MSN in striatum labeled with neurobiotin that we patched. (C) Representative traces obtained in striatal MSNs from *Trpv3*^{-/-} + Ctrl and *Trpv3*^{-/-} + TRPV3 mice obtained at RMP by whole-cell current-clamp recordings. (D) The summary of AP No. at RMP. AP No. of each current step was analyzed by two-tailed nonparametric Mann–Whitney test. (E) The statistics of RMP values and input resistance (IR) (F) values at RMP. (G) Representative traces of mEPSC in striatal MSNs from *Trpv3*^{-/-} + Ctrl and *Trpv3*^{-/-} + TRPV3 groups by whole-cell voltage-clamp recordings. (H) The frequency of mEPSC and associated cumulative probability (I). (J) Representative images showing the morphology of spines in MSNs from *Trpv3*^{-/-} + Ctrl and *Trpv3*^{-/-} + TRPV3 mice and the summary of spine density with a schematic of the dendrite. Exacerbation of neurological deficit scores (K) shown as the median with 95% CI and analyzed by Mann–Whitney test, brain water content (%) (L) and infarction volume (%) (M) in *Trpv3*^{-/-} + TRPV3 mice subjected to cerebral I/R (1.5/24 h) injury. Data are expressed as mean \pm SEM, n shows the number of neurons or the number of mice. * $P < 0.05$, ** $P < 0.01$ and *** $P < 0.001$. Also see Supporting Information Fig.

S4H for unnoticeable change of the complex dendritic trees in *Trpv3*^{-/-} + Ctrl and *Trpv3*^{-/-}+TRPV3 neurons and Table S5 for electrophysiological parameters of mouse striatal MSNs between *Trpv3*^{-/-} + Ctrl and *Trpv3*^{-/-} + TRPV3; and Table S6 for mEPSC parameters of mouse striatal MSNs between *Trpv3*^{-/-} + Ctrl and *Trpv3*^{-/-}+ TRPV3 groups.

Figure 6 Pharmacological inhibition of TRPV3 channels by natural compound forsythoside B reduces neural excitability and alleviates ischemia injury. (A) Raw traces from a single electrode recording of mouse primary cortical neurons in multi-electrode array (MEA) recordings. Cortical neurons produce fewer spikes after 100 $\mu\text{mol/L}$ FB treating. (B) Heat-maps of representative MEA recordings from cortical neurons culture at 37 °C. Application of FB decreases bust frequency of cortical neurons. (*n* shows the number of wells, * $P < 0.05$ and ** $P < 0.01$ versus before group). (C) Representative traces for mEPSCs in striatal MSNs from before and after 50 $\mu\text{mol/L}$ FB application in WT + Ctrl and WT + TRPV3 groups by whole-cell voltage-clamp recordings. (D) The frequencies of mEPSC with associated cumulative probability (E) before and after 50 $\mu\text{mol/L}$ FB application in WT + Ctrl and WT + TRPV3 groups (** $P < 0.01$ and *** $P < 0.001$, paired student *t*-test). (F) Application of FB (50 $\mu\text{mol/L}$) induced hyperpolarized RMP of striatal MSNs from WT + Ctrl and WT + TRPV3 groups. (G) The morphology of primary mouse cortical neurons before and after OGD or in the presence and absence of different concentrations of FB or edaravone. Scale, 100 μm . Summary of cell viability of primary mouse cortical neurons subject to OGD/R injury. *n* represents the number of independent experiments, * $P < 0.05$, ** $P < 0.01$, versus OGD group, one-way ANOVA with the Geisser–Greenhouse correction followed by uncorrected Fisher’s LSD test from GraphPad 8. (H) Representative images of TTC-stained mouse brain slices at 24 h after reperfusion. The white regions indicate the infarct size, and the red regions indicate the viable tissues. (I) Reduction of infarction volume (%) in mice subjected to cerebral I/R (1.5 h/24 h) injury by intravenous injection of FB (10, 20, and 40 mg/kg) at the beginning of reperfusion. (J) Decreases of neurological deficit scores in mice subjected to cerebral I/R injury by intravenous injection of FB (10, 20, and 40 mg/kg) or 3 mg/kg edaravone at the beginning of reperfusion. (K) Lack of neuroprotection by FB (10 mg/kg), neurological deficit scores (L) and brain water content (M) after I/R injury in *Trpv3* null mice by intravenous injection at the beginning of reperfusion. *n* indicates the number of mice,

* $P < 0.05$, ** $P < 0.01$ and *** $P < 0.001$ versus vehicle group. Data for infarct volume were analyzed by one-way ANOVA followed by Dunnett's test. Data for neurological deficit scores are shown as the median with 95% CI and analyzed by Kruskal–Wallis test followed by Dunn's test from GraphPad 8. Other data are expressed as mean \pm SEM. Also see Supporting Information Fig. S5 for pharmacological inhibition of TRPV3 channels by forsythoside B (FB) that reduces neural excitability, and Table S7 for effects of FB (50 $\mu\text{mol/L}$) on electrophysiological parameters and Table S8 for effects of FB (50 $\mu\text{mol/L}$) on mEPSC parameters of mouse striatal MSNs from WT + Ctrl and WT + TRPV3 groups.

Figure 7 A proposed mechanism for a causative role of overactive intracellular proton-sensitive and Ca^{2+} -permeable TRPV3 channels in brain ischemic injury. Upon cerebral ischemia injury occurs, quick energy depletion leads to metabolic acidosis resulted from metabolizing pyruvate into lactic acid through the anaerobic pathway. Accumulation of lactic acid^{1,90} results in intracellular rise of H^+ that activates TRPV3 channels likely through binding to intracellular N-terminal H426 site for induction of calcium influx²⁴. Intracellular calcium also activates TRPV3 channels through binding to R696 site⁹¹. Energy depletion induces anoxic depolarization that can activate some voltage-dependent calcium channels. In addition, the extracellular acidification ($\text{pH}_o = 6.5$)⁹² can activate acid-sensing ion channels (ASICs) and potentiate the activation of transient receptor potential melastatin 7 (TRPM7) channels for further Ca^{2+} toxicity²³. All these cascade events including calcium releases from endoplasmic reticulum and mitochondria cause intracellular calcium overload, thus leading to neuronal death. Conversely, silencing of pharmacological inhibition of overactive TRPV3 channels by forsythoside B reduces neuronal death and alleviates ischemic brain injury.

