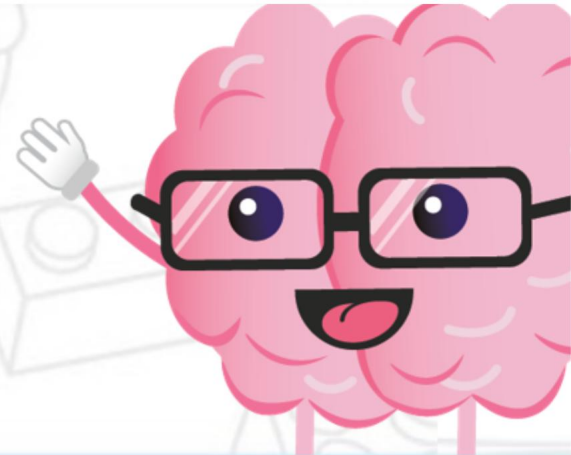
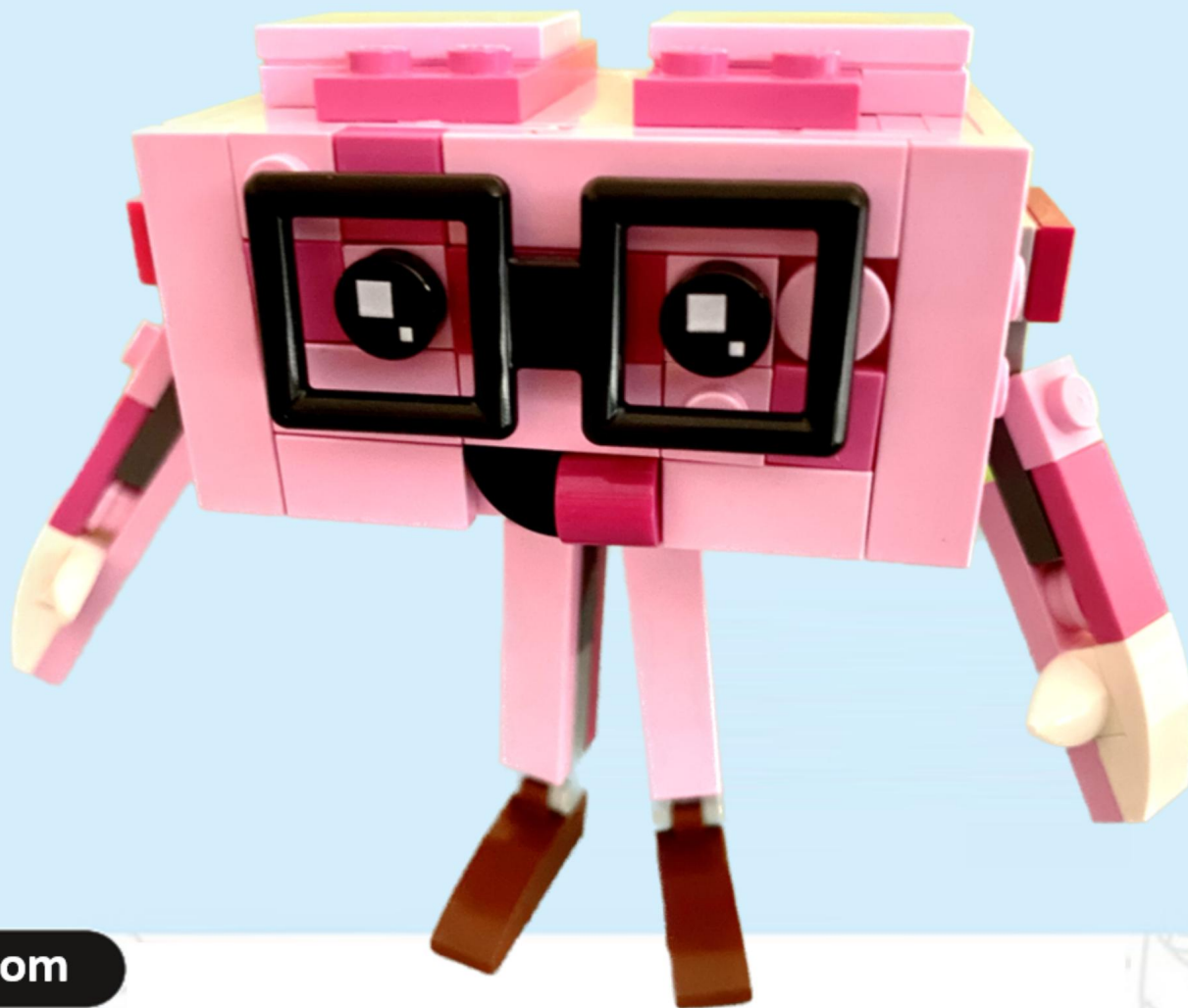




**BIO  
FRIENDS™**



**96768  
Mini-Brain**





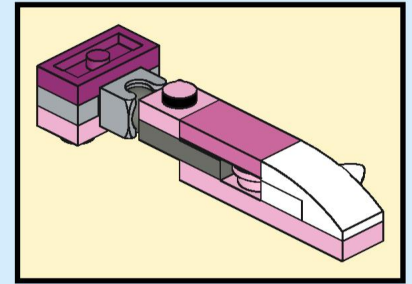
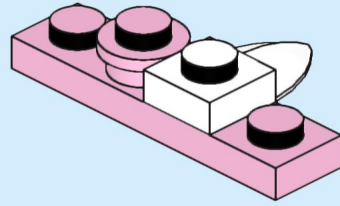
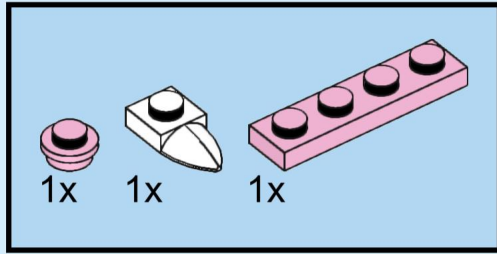
# Making **organoid cell** models can be challenging

Axion BioSystems offers easy-to-use tools that enable consistent organoid production and provide deep insights into their biology.

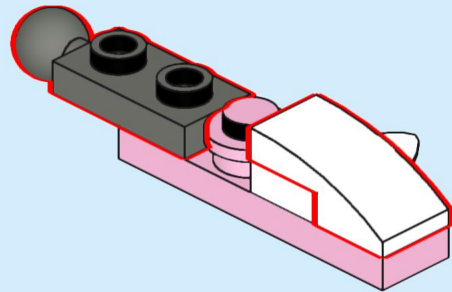
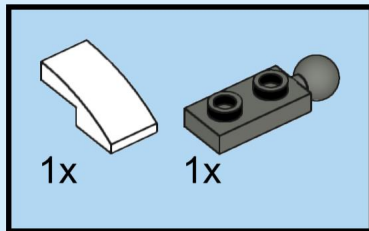
Excited to learn more about the future of organoid-based research?  
**Scan the QR code.**



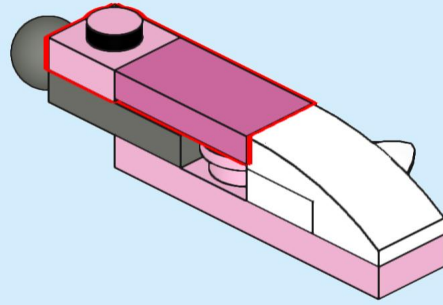
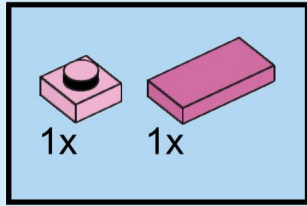
# 1



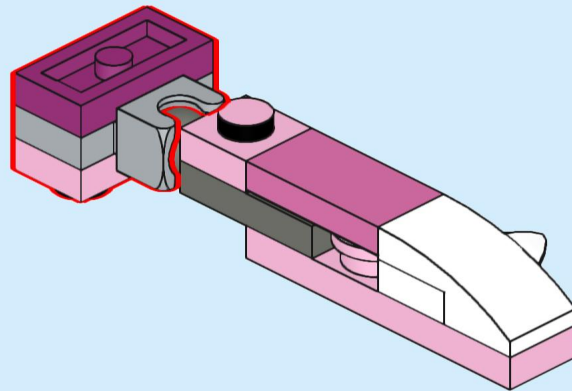
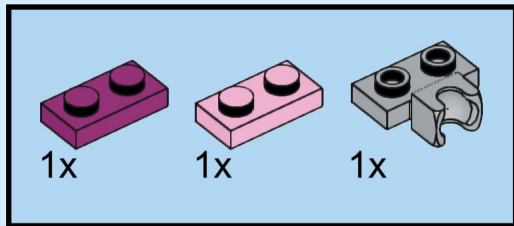
# 2



3

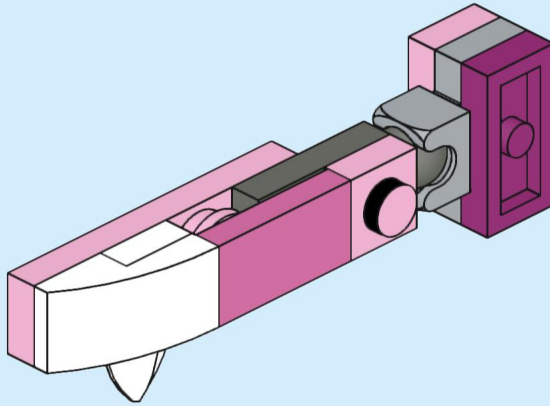


4

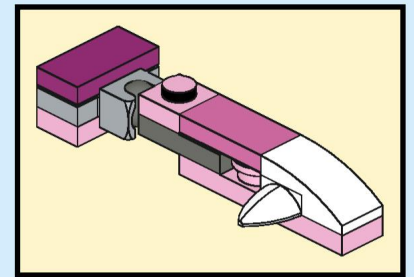
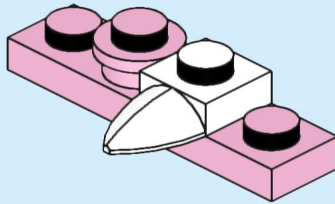
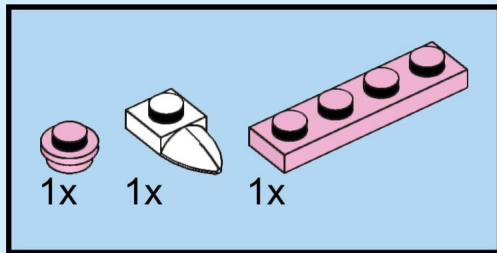




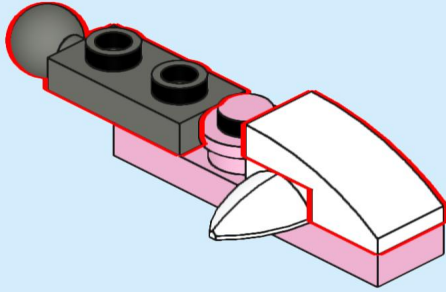
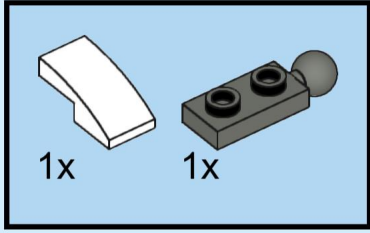
# 5



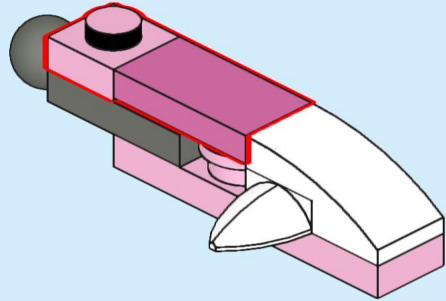
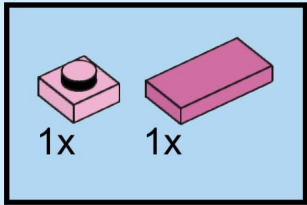
# 6



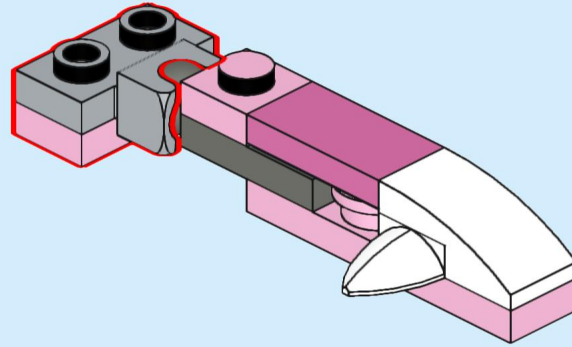
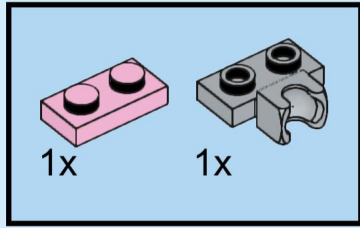
7



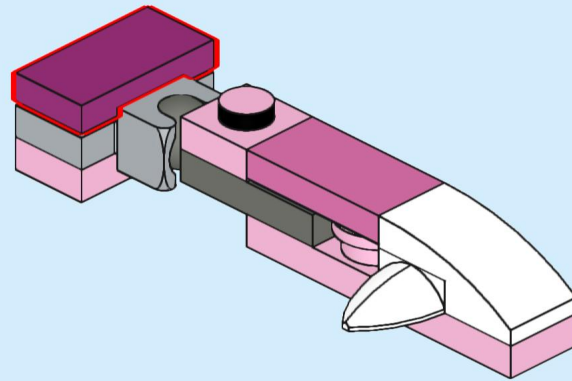
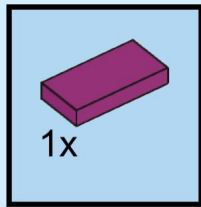
8



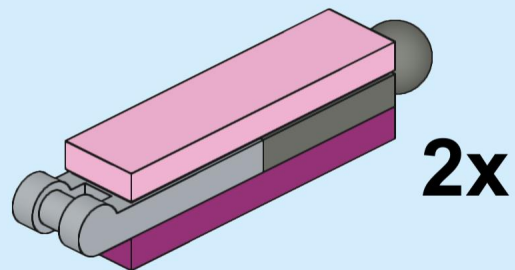
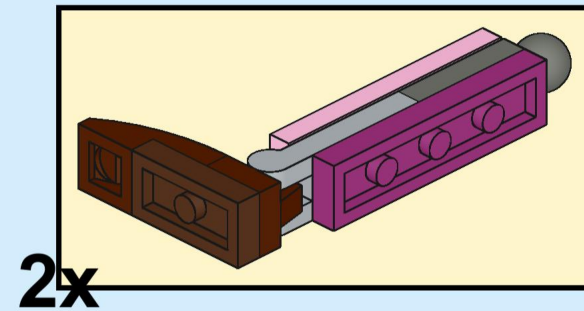
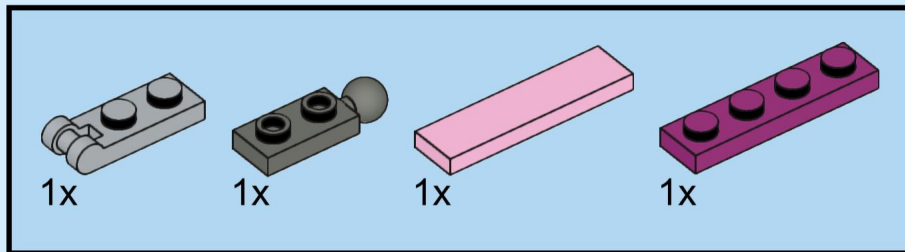
# 9



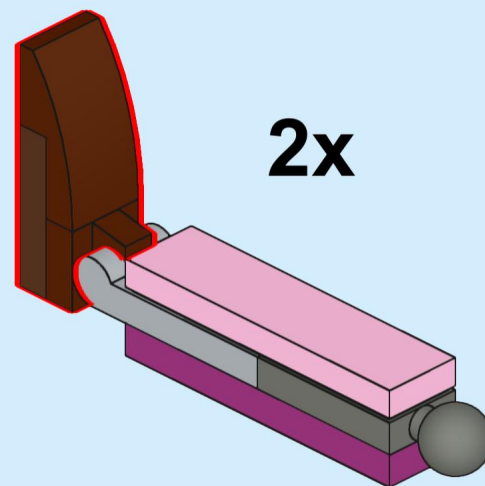
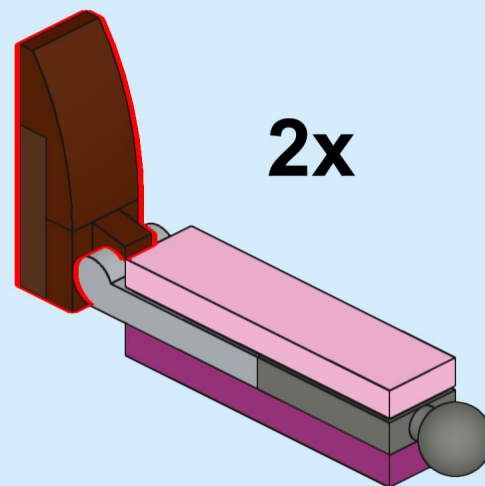
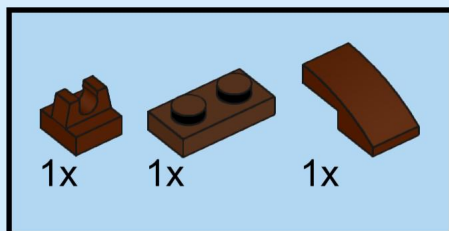
# 10



# 11

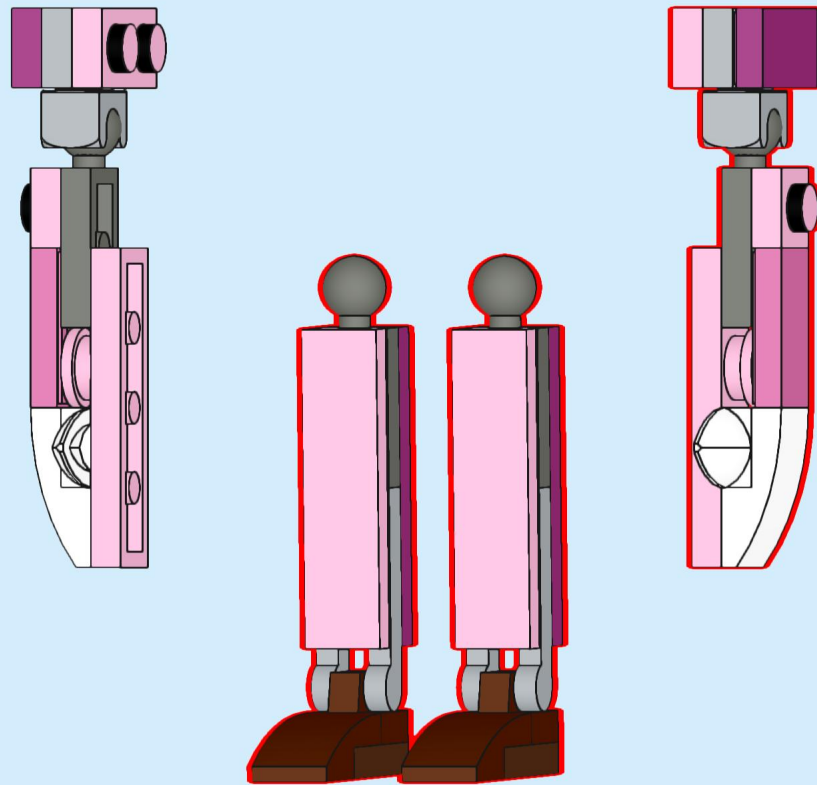


# 12

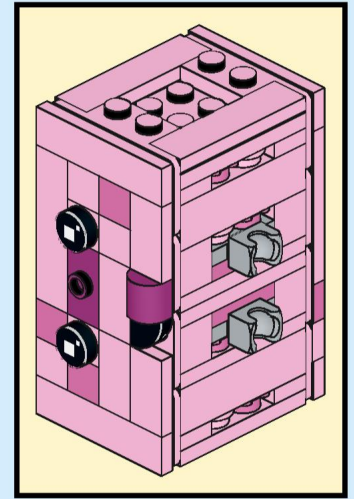
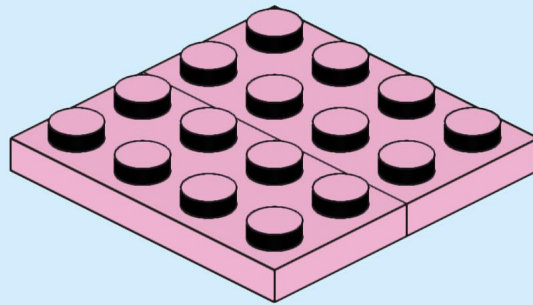
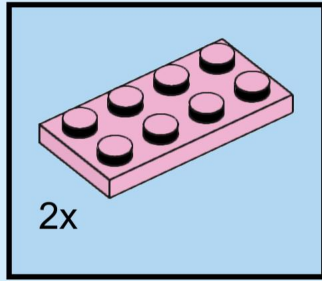




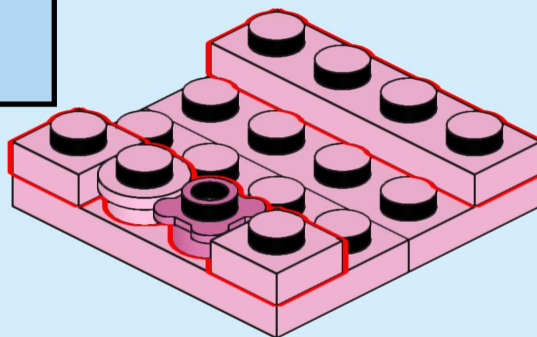
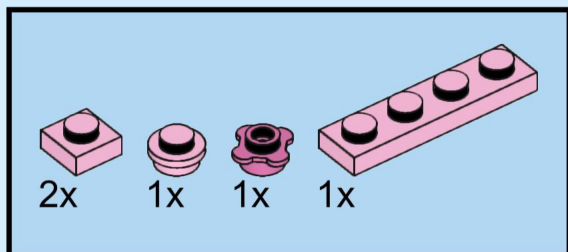
13



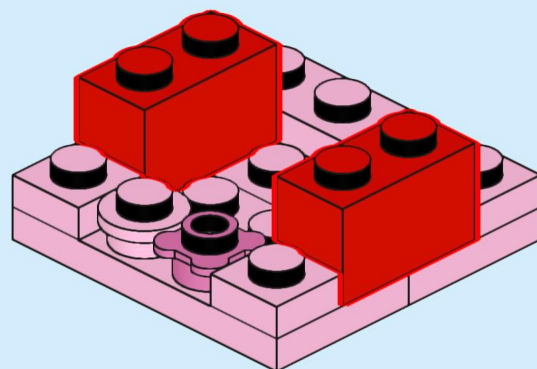
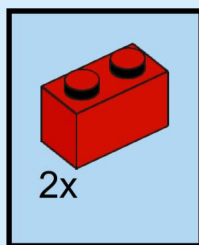
# 14



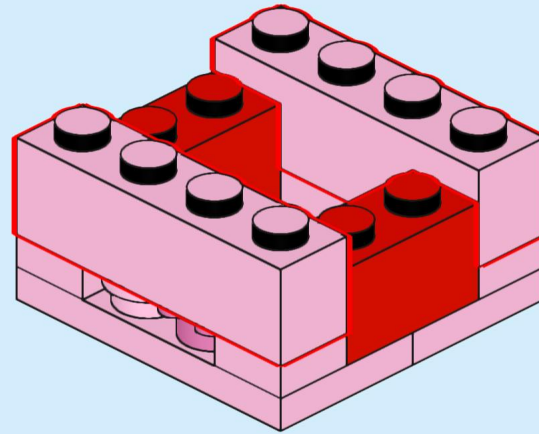
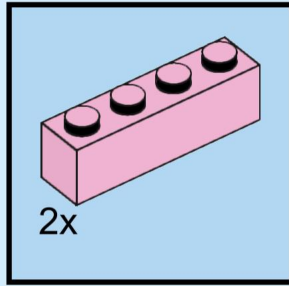
# 15



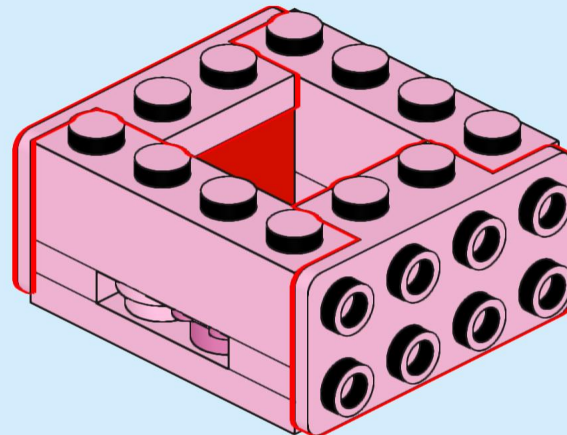
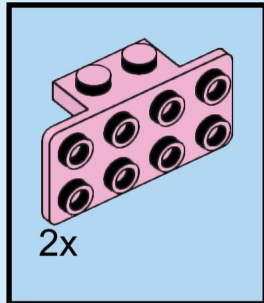
# 16



# 17

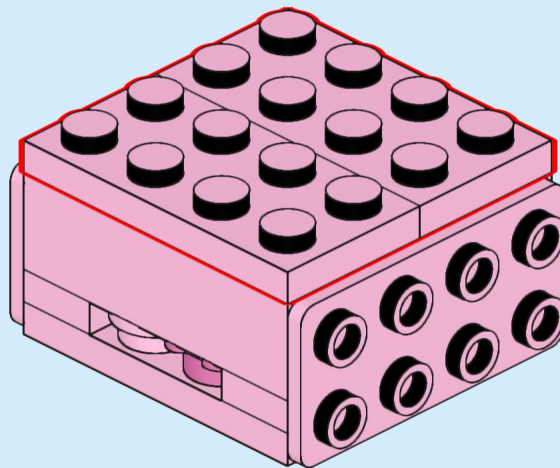
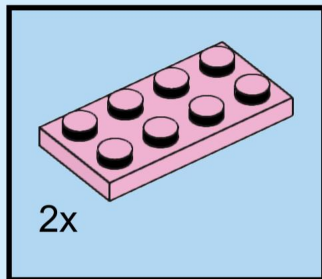


# 18

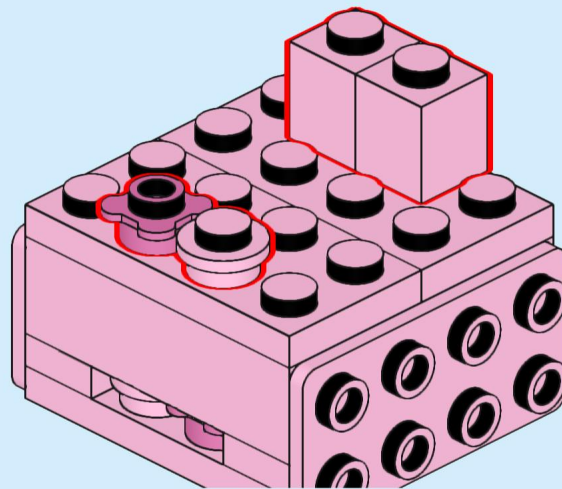
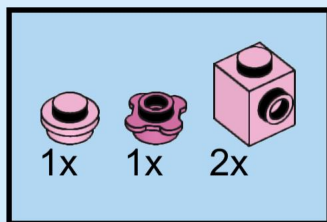




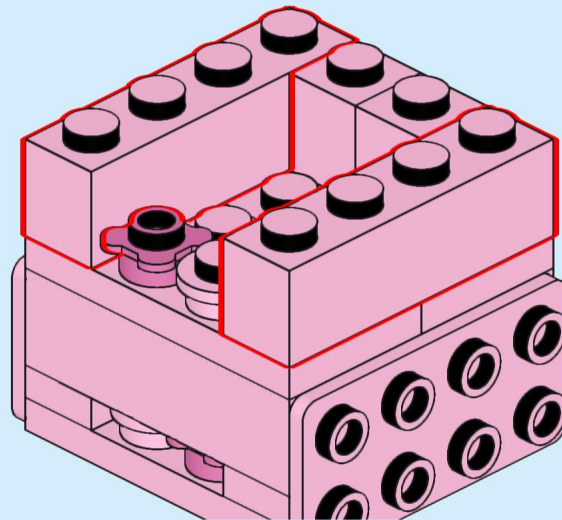
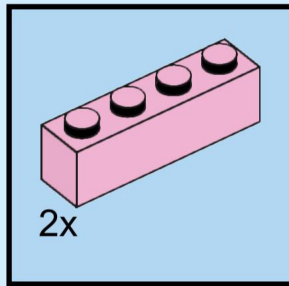
# 19



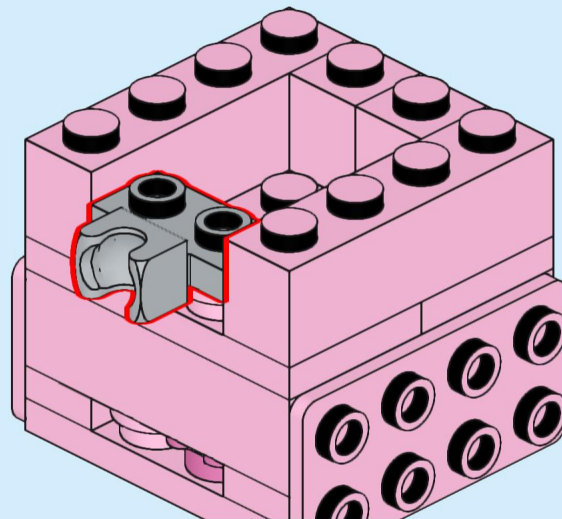
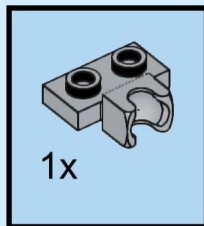
# 20



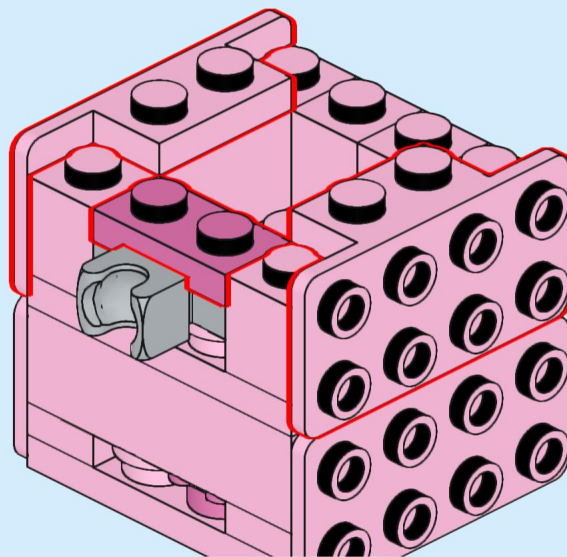
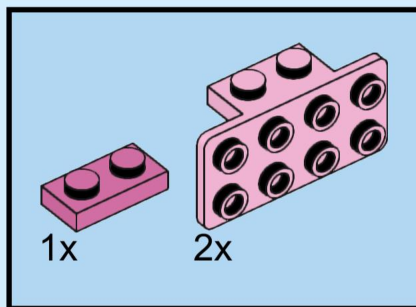
# 21



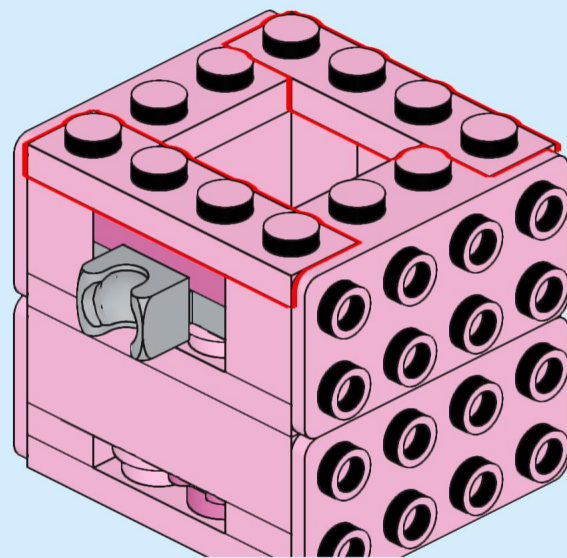
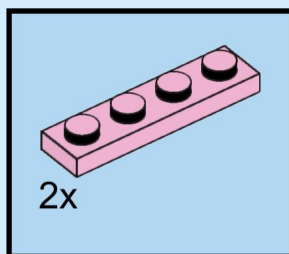
# 22



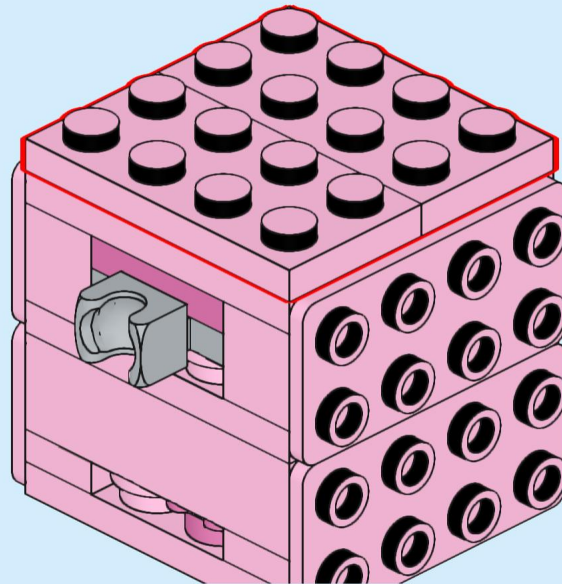
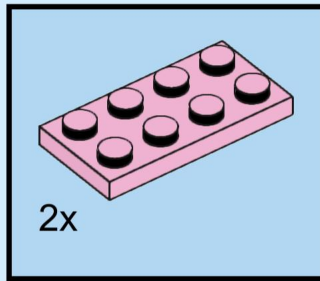
# 23



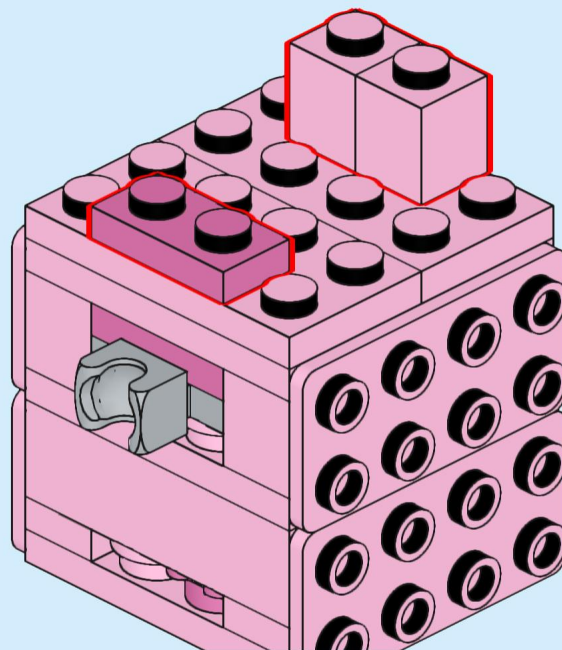
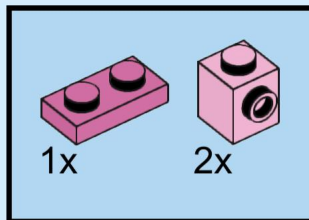
# 24



# 25

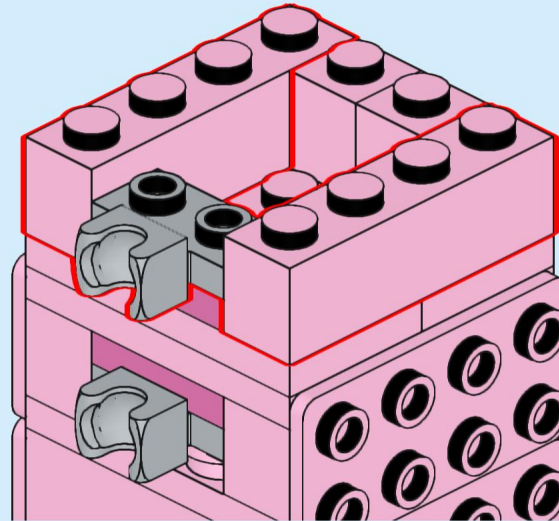
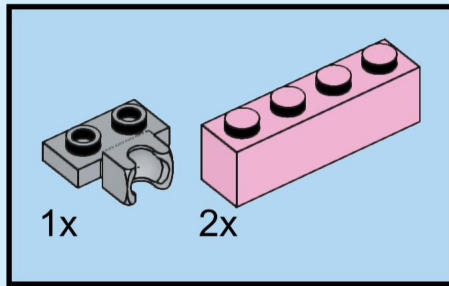


# 26

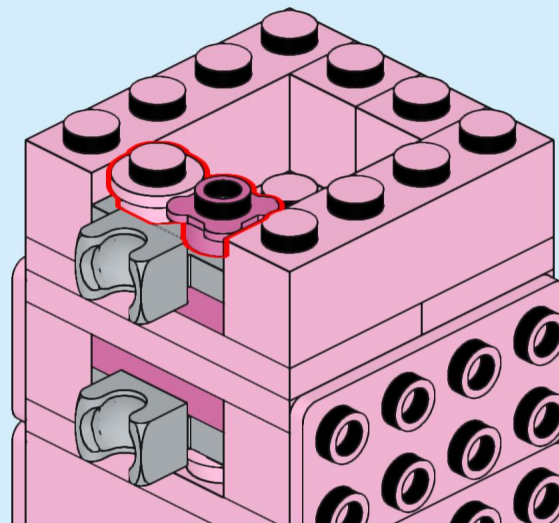
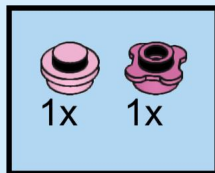




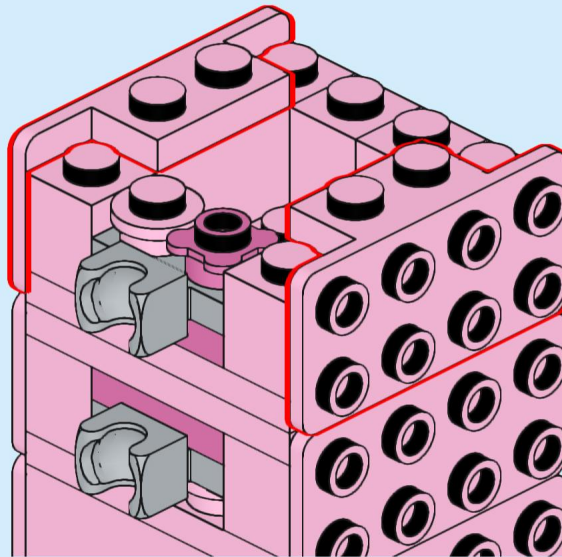
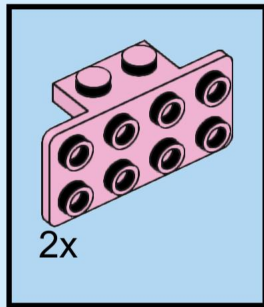
# 27



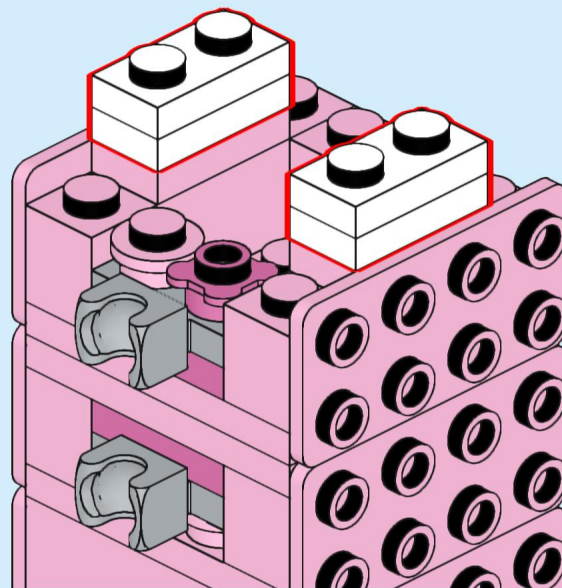
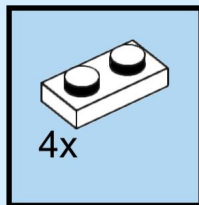
# 28



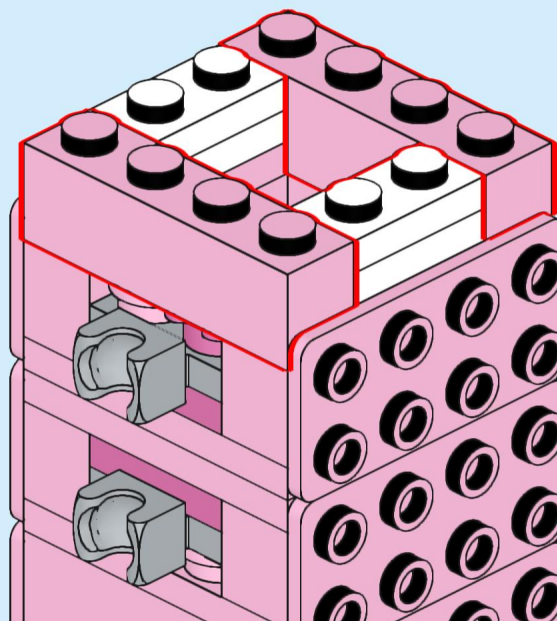
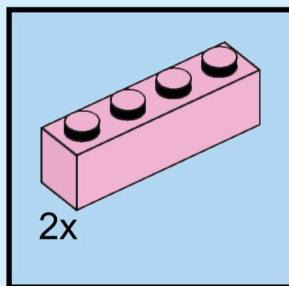
# 29



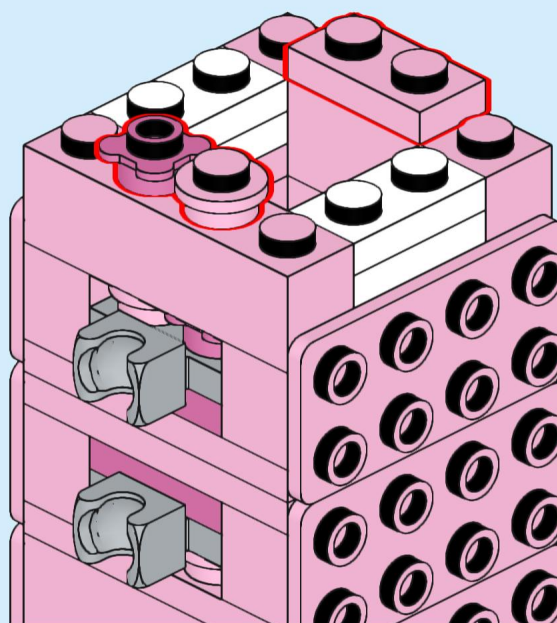
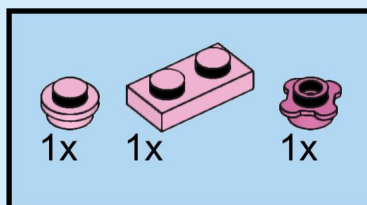
# 30



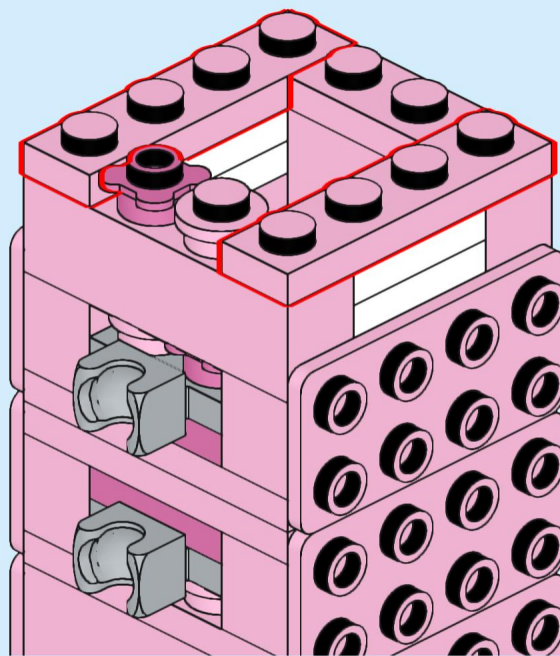
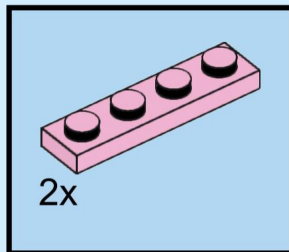
# 31



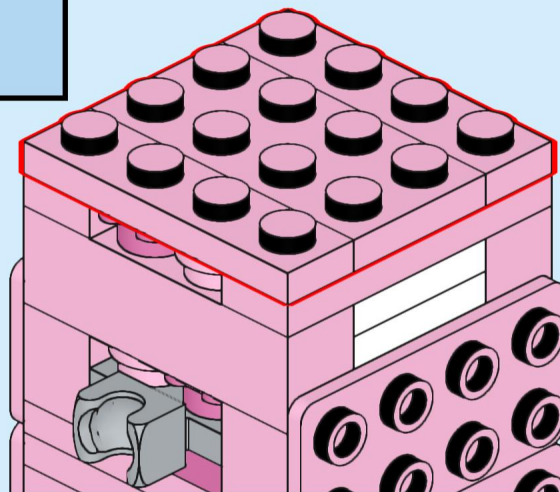
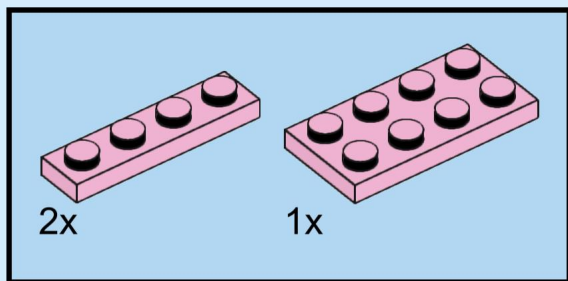
# 32



# 33

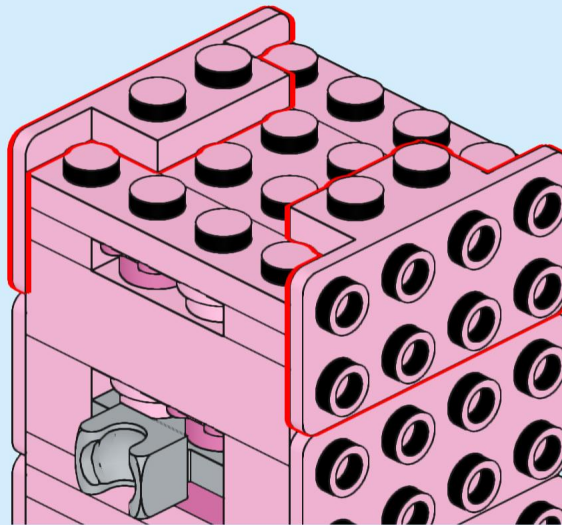
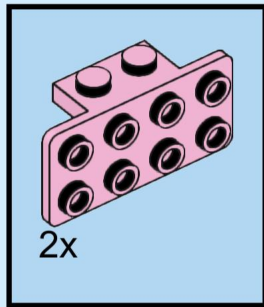


# 34

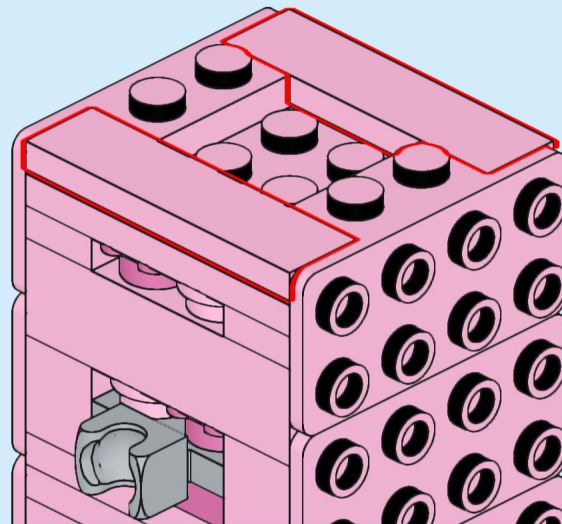
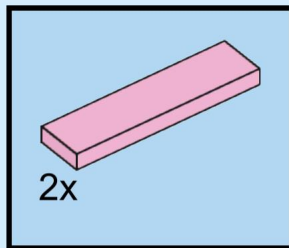




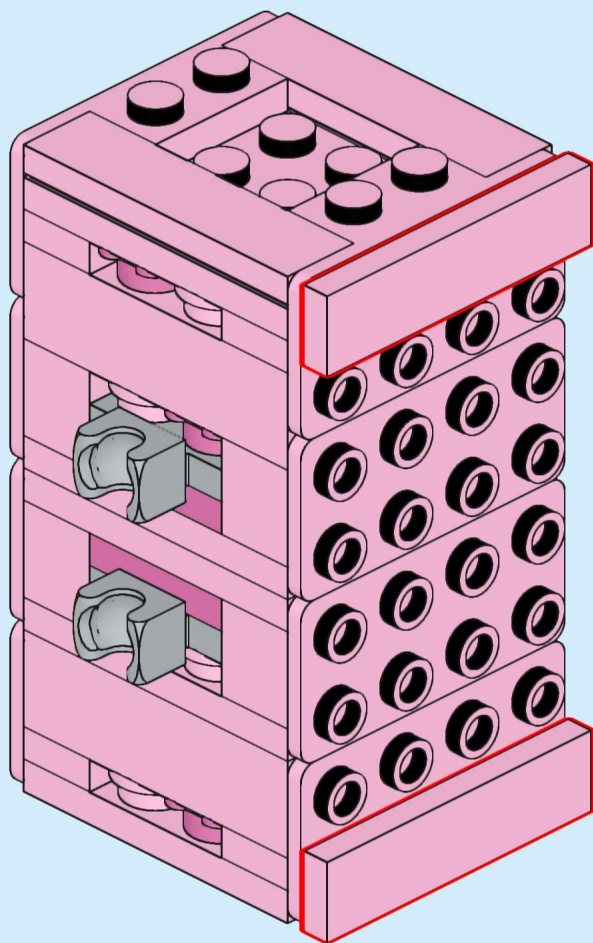
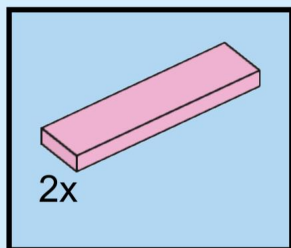
# 35



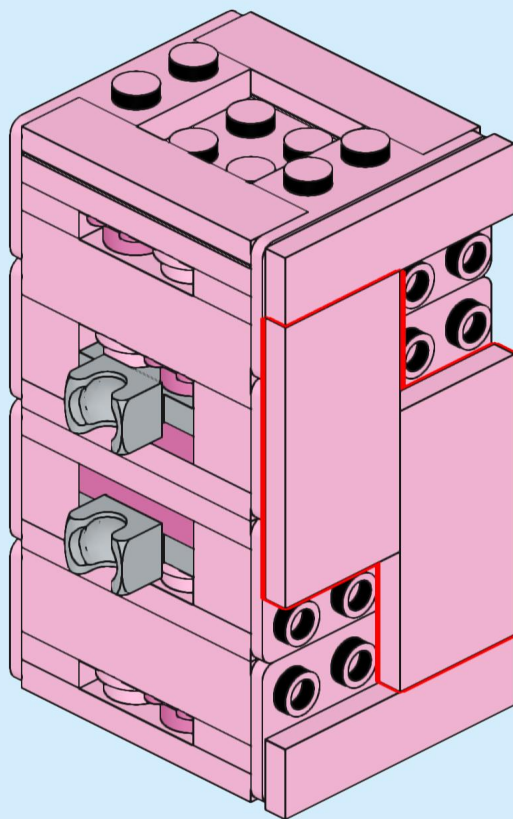
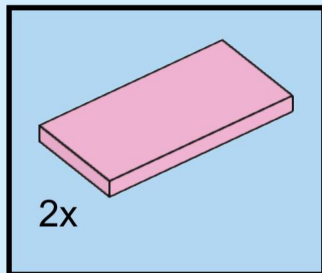
# 36



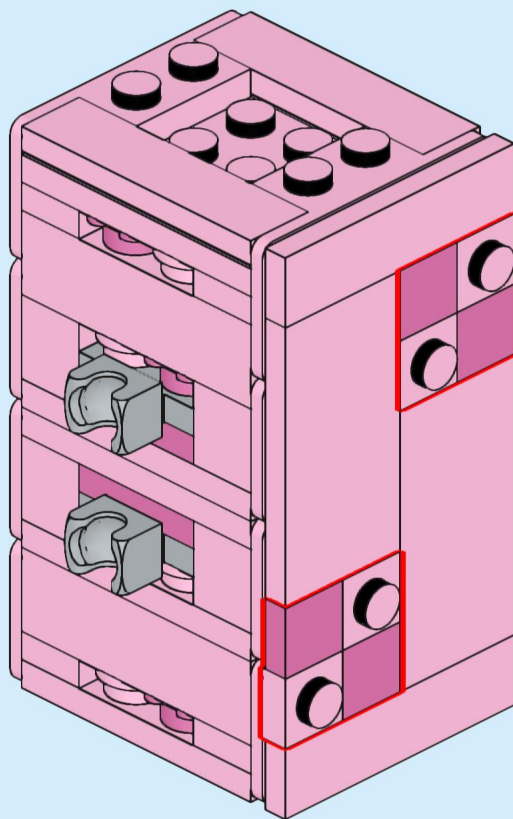
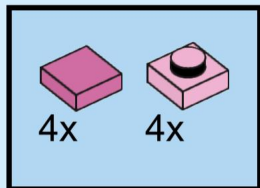
# 37



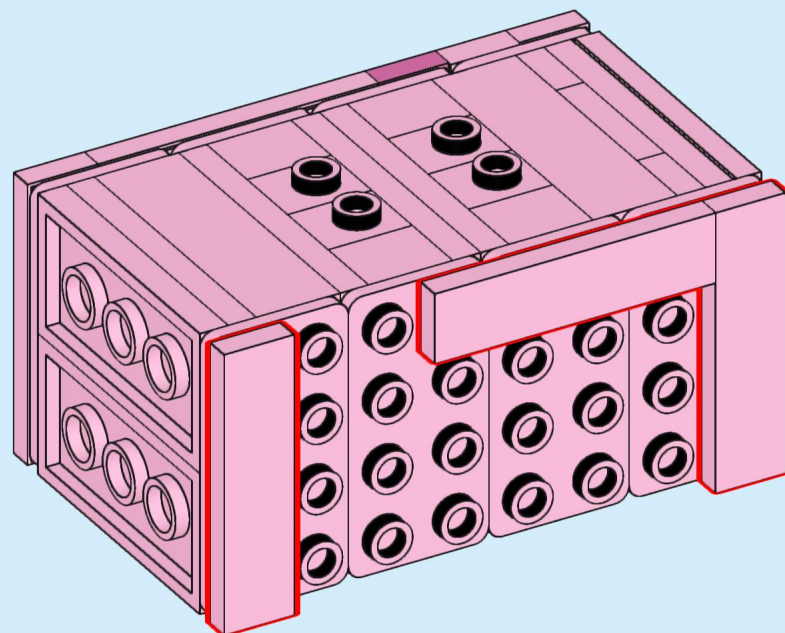
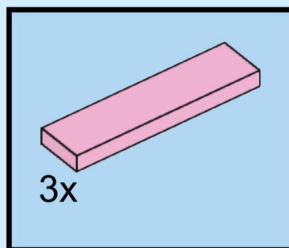
# 38



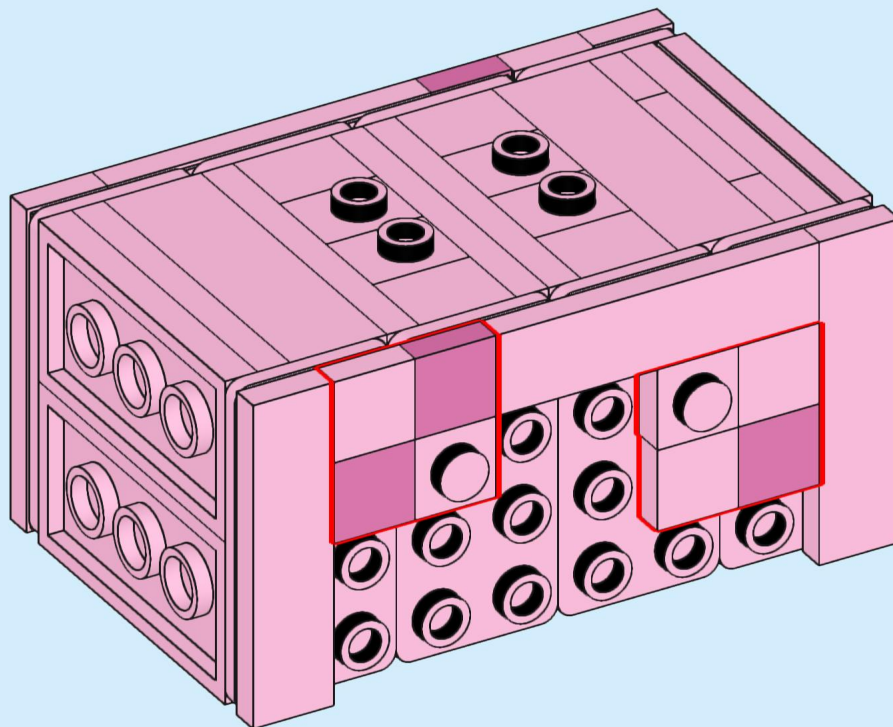
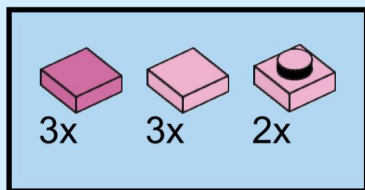
# 39



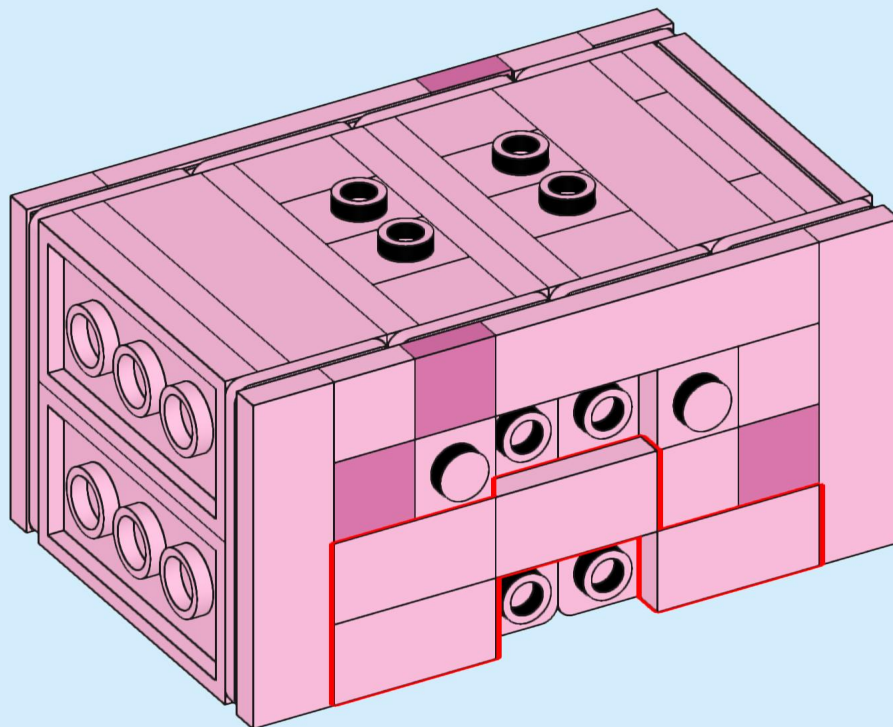
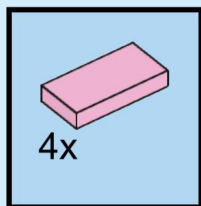
40



# 41

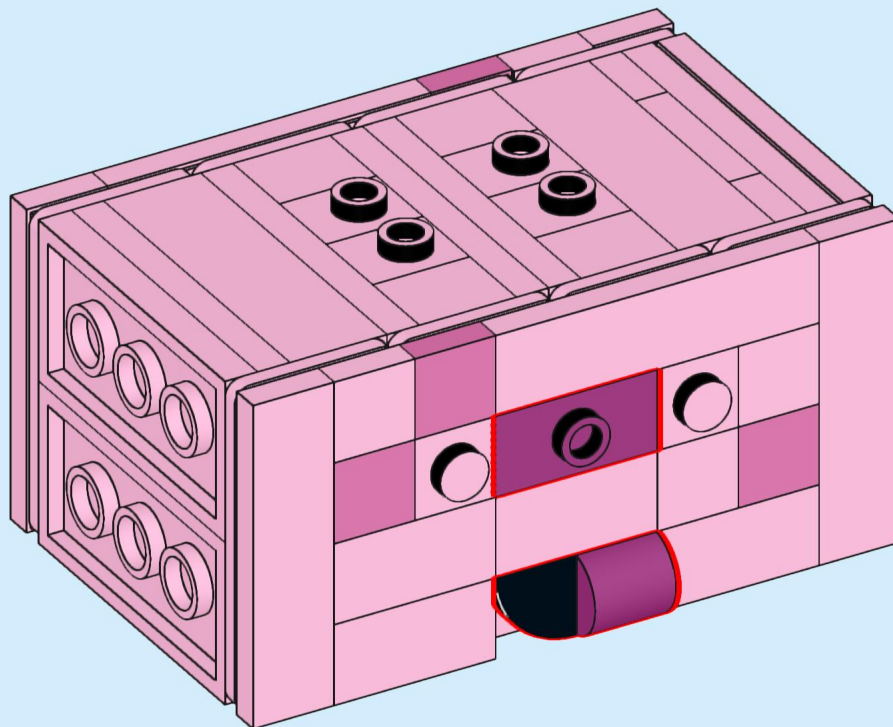
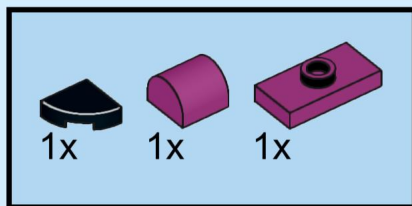


# 42

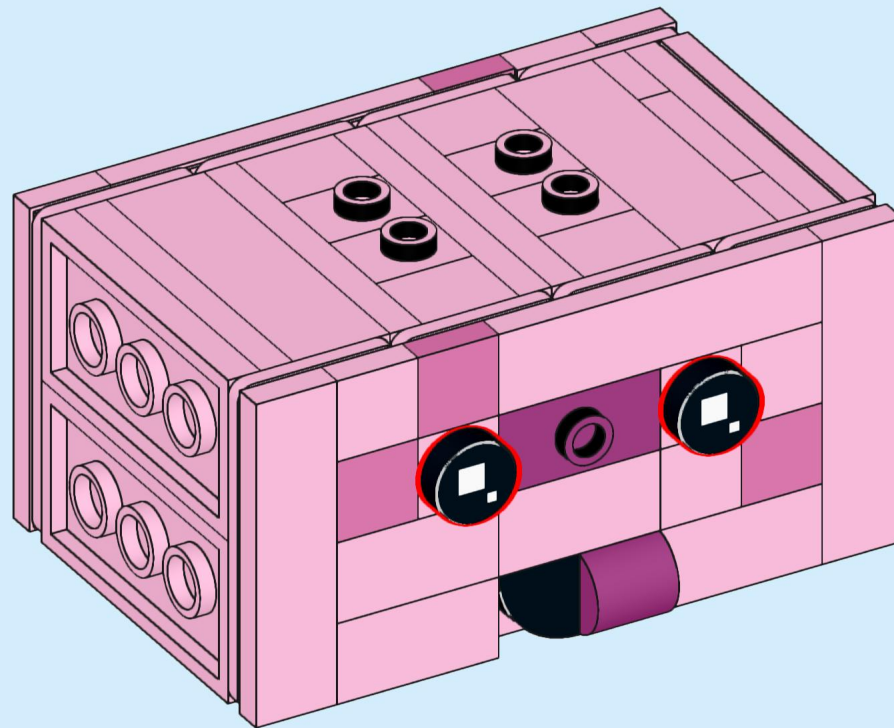




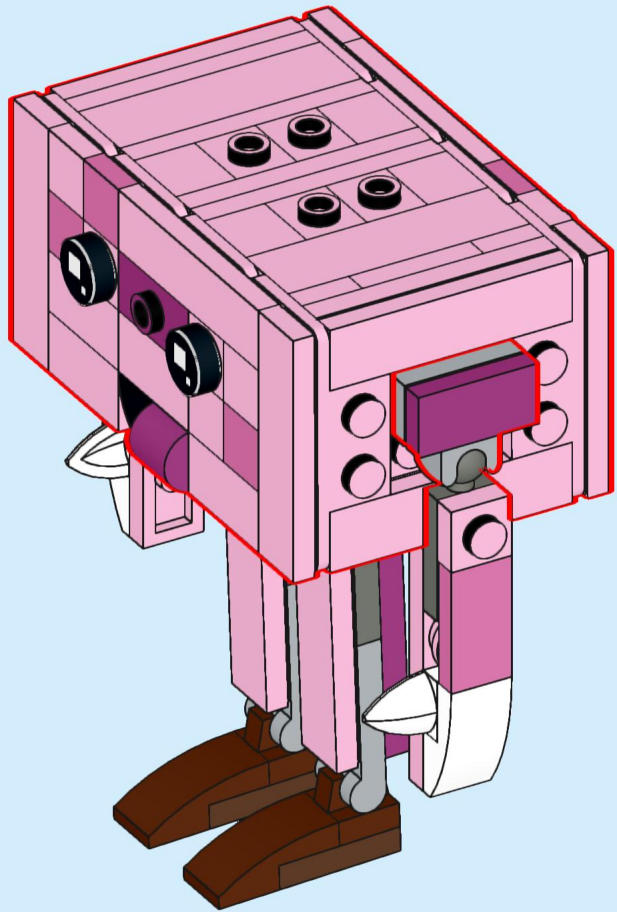
# 43



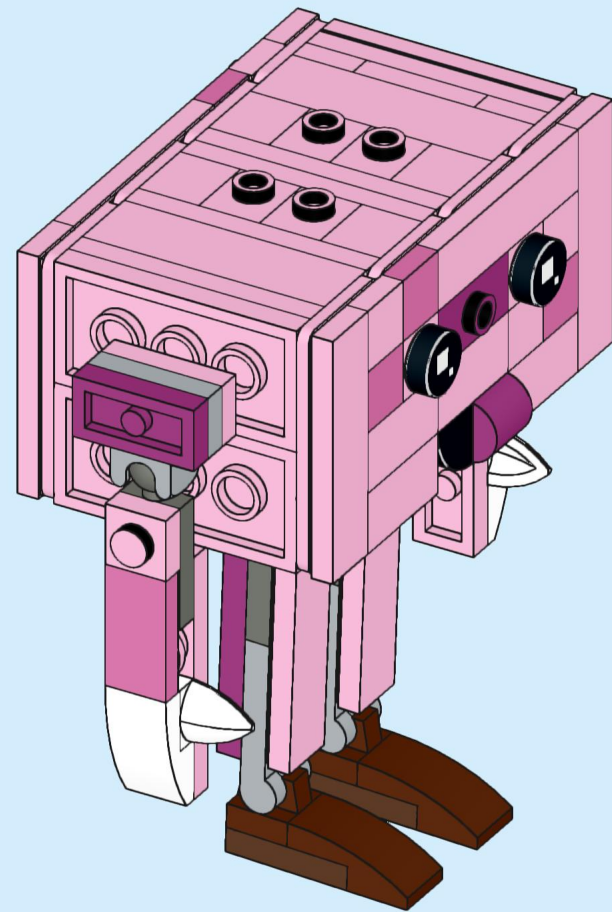
44



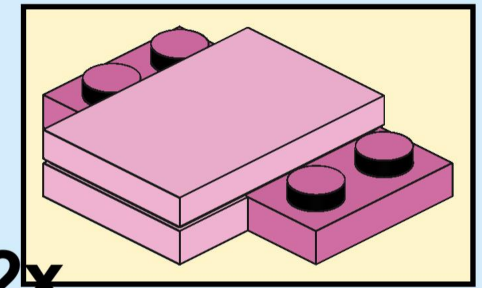
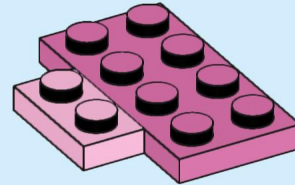
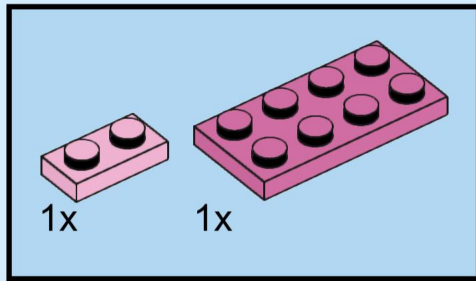
45



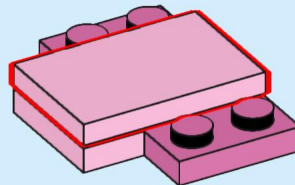
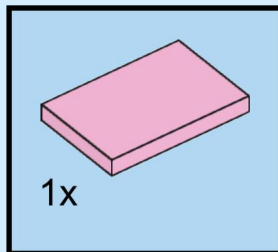
46



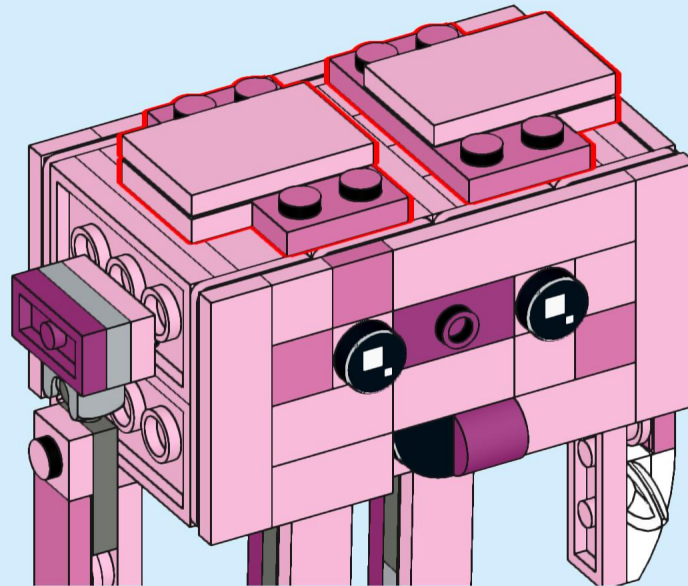
# 47



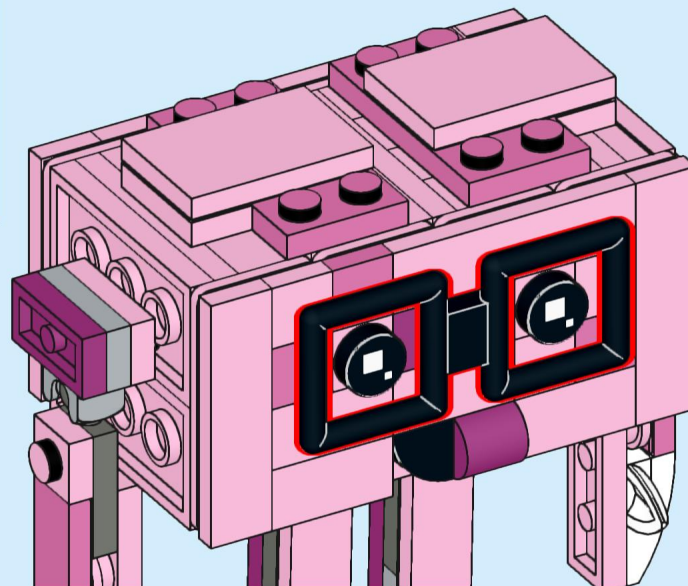
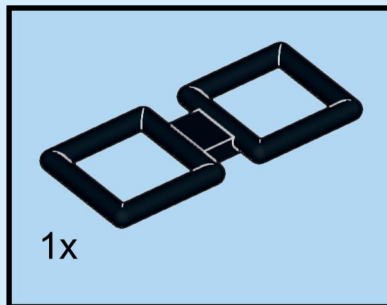
# 48



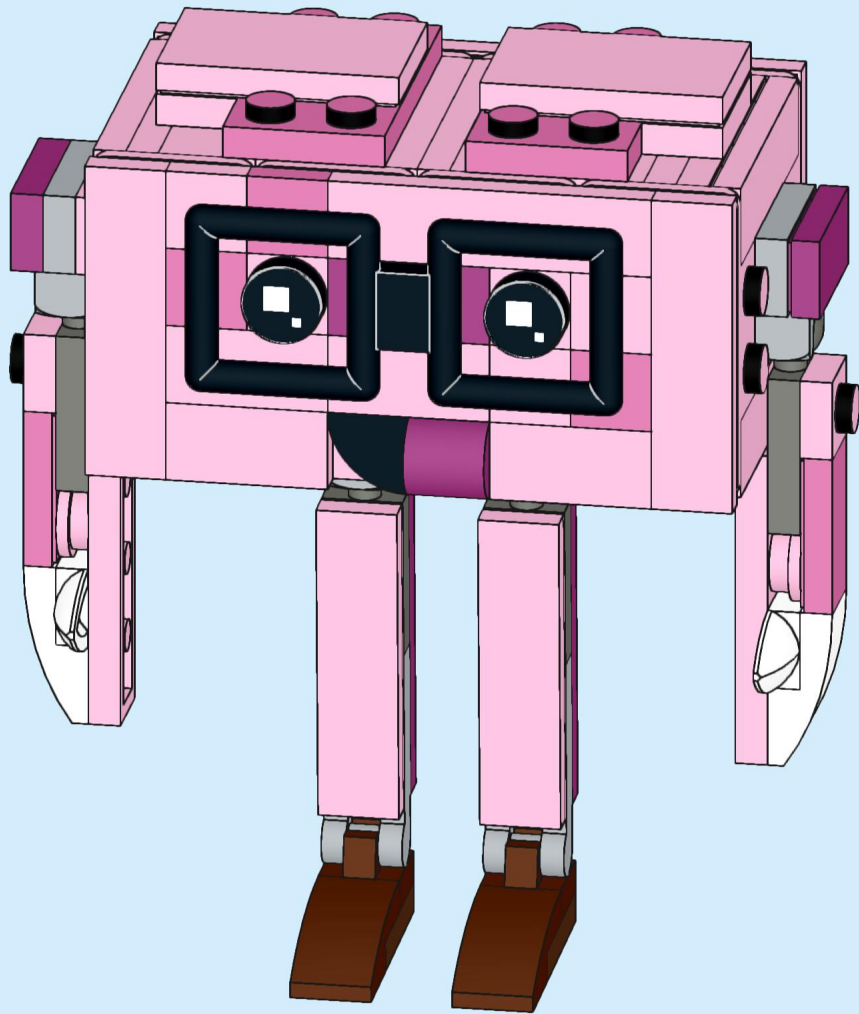
# 49



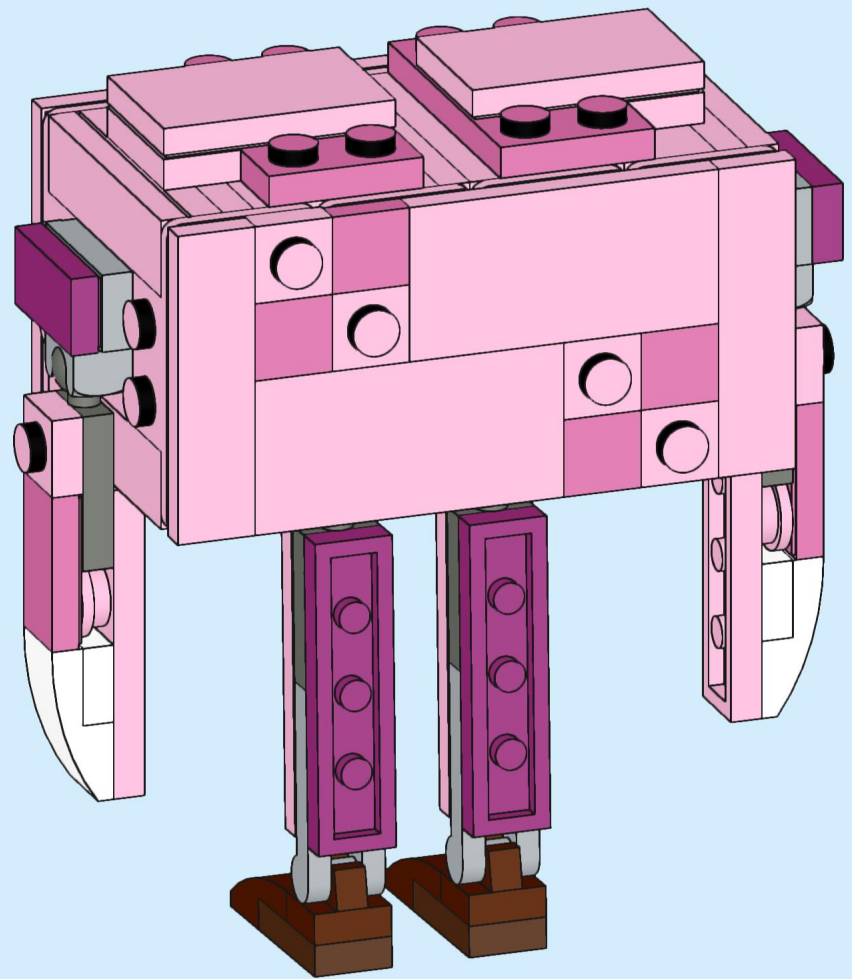
# 50



# 51



# 52



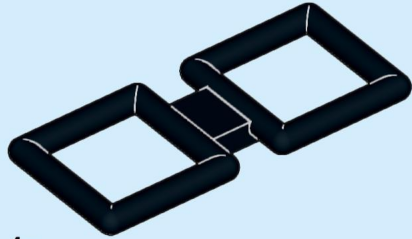




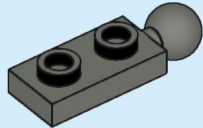
2x  
6284077



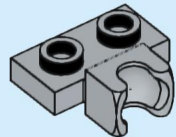
1x  
6172383



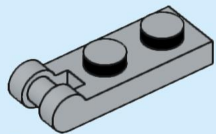
1x  
6208551



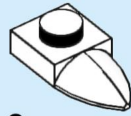
4x  
6123814



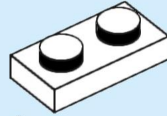
4x  
6043656



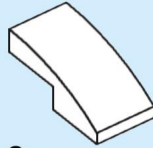
2x  
4515369



2x  
4224792



4x  
302301



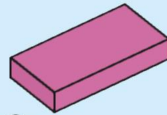
2x  
6034044



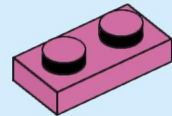
7x  
6133726



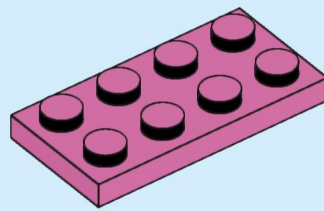
4x  
6170298



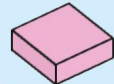
2x  
6056381



2x  
6057387



2x  
6056263



3x  
6251940



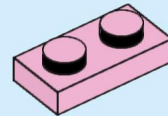
6x  
4517996



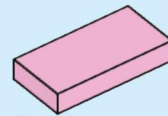
10x  
6031883



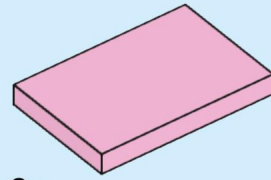
4x  
4621554



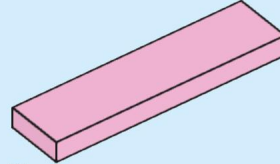
5x  
4654128



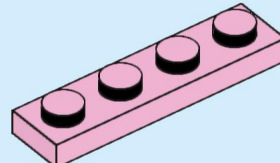
4x  
4580010



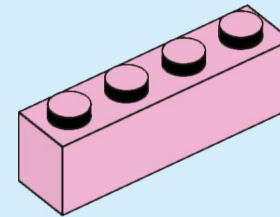
2x  
6290982



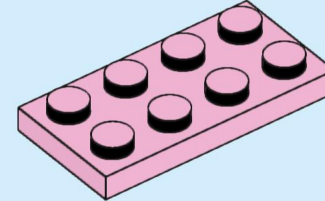
9x  
4621552



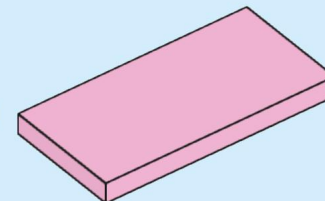
9x  
6002148



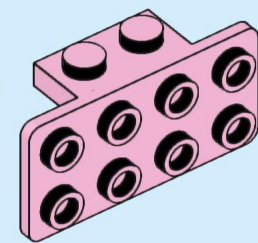
8x  
4518890



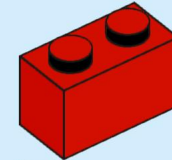
7x  
6204535



2x  
6106713



8x  
6122589



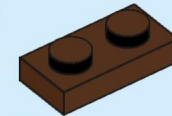
2x  
300421



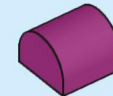
2x  
6030720



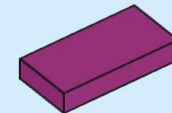
2x  
6069984



2x  
302325

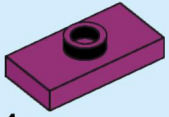


1x  
6278442

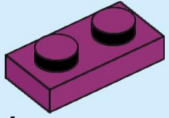


1x  
4611415

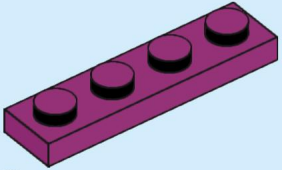




1x  
6101073



1x  
6103415



2x  
6037652





For research use only. Not for use in diagnostic procedures.