

Drug Dosing Calculator



Contents

Chapter 1. Overview	3
1.1. Drug Dosing Calculator Overview	3
1.2. Stock Solution Conditions	3
1.3. Target Dose Conditions.....	3
1.4. Well Conditions	4
1.5. Solution Results	4
Chapter 2. Definitions.....	5
2.1. Starting Conditions.....	5
2.2. Target Conditions	5
2.3. Well Conditions	5
2.4. Results	6
Chapter 3. Equations.....	7
3.1. Solution Design	7



Chapter 1. Overview

1.1. Drug Dosing Calculator Overview

Axion's Drug Dosing Calculator calculates the concentrations and volumes needed to prepare dosing solutions for plate-based experiments.

1.2. Stock Solution Conditions

1. Enter your concentrated stock solution (**Stock**) and the vehicle concentration in the stock solution (**Vehicle**).
 - 1.1 Ensure proper units are selected using the dropdown option.
 - 1.2 *For example, a compound prepared in all DMSO would typically need to be entered as 100% DMSO.*

Step 1

Starting Stock Conditions:

[Stock] *	Units	Vehicle	[Vehicle] *
	mM	DMSO	100 %

1.3. Target Dose Conditions

1. Enter your target cell plating density in the **Cells per Well** field, as well as the cell plating volume (**V_{plating}**) and number of wells to be plated.
2. Multiple plating densities may be added or removed with the **plus** and **minus** buttons.
 - 2.1 **Note:** Order the list of dosage concentrations with proper units from highest to lowest.
 - 2.2 Additionally, Spare Volume provides room to account for extra solution in each aliquot due to pipetting loss or dead volume.

Step 2

Target Dose Conditions:

Group	[Dose]	Units	Replicates	Remove
1		μM	1	

+ Add Dose Row

Spare Volume

%



1.4. Well Conditions

1. Enter your well-starting solution volume (**Starting Well Volume**) and the % of desired final vehicle concentration into **Final [Vehicle]**.

1.1. **Dose Volume** represents the volume of drug solution added to the well.

2. To add to the initial volume or replace an equivalent amount of volume with dose solution, click on **Add To** or **Replace** respectively.

Step 3

Experiment Setup:

Well Contents sodium	Starting Well Volume 100 μL	Final [Vehicle] 0.1 %
Dose Volume 10 μL	Dose Method ☒ Add to ☐ Replace Add to: The dose solution is added to the existing well volume. Replace: The same volume is removed from the well and replaced with dose solution.	

1.5. Solution Results

Then, click **Calculate Results**. The following outputs are provided:

1. **Product** identifies the solution being prepared.
2. **[Conc]** provides the target concentration of the product.
3. Solution identifies the source solution used to prepare the product.
4. **V_{solution}** provides the required volume of the source solution.
 - 4.1. It is rounded to the nearest 0.1 μL and is not allowed to be less than 10 μL . V_{diluent} is adjusted accordingly.
5. **Diluent** identifies the diluent used to prepare the product.
6. **V_{diluent}** provides the required volume of diluent.

For each solution (Product) you need to make, add **V_{diluent}** to **V_{solution}** slowly, mixing as needed.

Note: Adding **V_{solution}** to **V_{diluent}** may cause solubility issues. Always dilute high to low.

Note: The starting stock or vehicle concentration is considered too low if the desired final vehicle concentration cannot be achieved.

▼ Working Solution Preparation

Product	[Conc]	V _{solution}	Solution	V _{diluent}	Diluent
Work 1	55 μM	10 μL	Int 1	80.9 μL	Dil 1



Definitions

2.1. Starting Conditions

Stock Solution Conditions

[Stock]	Concentration of the stock solution
Vehicle	Vehicle control
[Vehicle]	Concentration of the vehicle in the stock solution

2.2. Target Conditions

Target Dose Conditions

[Dose]	Final concentration in the well
Replicates	Number of wells to receive the dose
Spare Volume	Percentage of overfill each aliquot will get to avoid running short.

2.3. Well Conditions

Experiment Setup

Well Contents	Solution or medium present in the well before dosing.
Starting Well Volume	Volume of solution or medium present in each well before dosing.
[Vehicle]	Final concentration of the vehicle in percent
Dose Volume	Volume of drug solution added to the well
Dose Method	Method of dose addition



2.4. Results

Results

Product	The final solution being prepared
[Conc]	Final concentration of the solution being prepared
V_{solution}	Required volume of the source solution
Solution	Source solution used to prepare the product
V_{diluent}	Required volume of diluent
Diluent	Diluent used to prepare the product



Chapter 3. Equations

3.1. Solution Design

Solution Design Calculations:

(n = number of drug solutions or length of the array)

$$\text{Number of Intermediates} = \text{Round} \frac{(\text{Log}(\text{Concentration}_{\text{stock}}) - \text{Log}(\text{Concentration}_{\text{work}_1}))}{2}$$

$$\text{Concentration}_{\text{working}_1} = \frac{(\text{Concentration}_{\text{dose}_n} * V_{\text{well}_{\text{final}}})}{V_{\text{add}}}$$

$$\text{For Add: } V_{\text{well}_{\text{final}}} = V_{\text{well}} + V_{\text{add}}$$

$$\text{For Replace: } V_{\text{well}_{\text{final}}} = V_{\text{well}}$$

$$\text{If } \left(\frac{\text{Log}(\text{Concentration}_{\text{stock}}) - \text{Log}(\text{Concentration}_{\text{work}_1})}{2} - (x - 1) \geq 1 \text{ then } C_{\text{drug}_n} = \frac{C_{\text{drug}_{n-1}}}{100} \right)$$

$$\text{Else if } \frac{\text{Log}(\text{Concentration}_{\text{stock}}) - \text{Log}(\text{Concentration}_{\text{work}_1})}{2} - (x - 1) \geq 0.5 \text{ and } \leq 1 \text{ then } C_{\text{drug}_n} = \frac{C_{\text{drug}_{n-1}}}{10}$$

Volume Calculations

$$\text{For working solutions: } (V_{\text{add}} * \text{Rep} + V_{S_{n+1}}) * \left(1 + \frac{V_{\text{extra}}}{100}\right)$$

For intermediates:

$$\text{If } V_{f_n} \leq (1000 \text{ ul}) \text{ then } V_{f_n} = 1000$$

$$\text{Else } V_{f_n} = (V_{\text{add}} * \text{Rep} + V_{S_{n+1}}) * \left(1 + \left(\frac{V_{\text{extra}}}{100}\right)\right)$$

$$V_{S_n} = \frac{(C_{\text{drug}_n} * V_{f_n})}{(C_{\text{drug}_{n-1}})}$$

$$V_{d_n} = V_{f_n} - V_{S_n}$$

**Vehicle & Diluent Preparation Calculations:**

$$Concentration_{vehicle_{working}} = \frac{Concentration_{vehicle_{final}} * V_{well_{final}}}{V_{add}}$$

$$C_{diluent_1} = C_{vehicle_{working}} - C_{vehicle_n} \text{ where } n = \text{the first solution where } C_{vehicle_n} \leq C_{vehicle_{working}}$$

$$Vf_{diluent_1} = Vd_n * \left(1 + \left(\frac{V_{extra}}{100}\right)\right) \text{ where } n = \text{the first solution where } C_{vehicle_n} \leq C_{vehicle_{working}}$$

$$VS_{diluent_1} = \frac{C_{diluent_1} * Vf_{diluent_1}}{100}$$

$$Vd_{diluent_2} = Vf_{diluent_2} - VS_{diluent_2}$$