

How to Use the Cell & Reagent Calculator for the Nucleofector™ 96-well Shuttle™ Technical Information

www.lonza.com/cell-reagent-calculator

1) Fill in the Top 2 Sections Based on Your Experimental Set Up and Stock Solutions

a) General

- Fill in the number of transfections you wish to perform, i.e. how many wells you will use. It is generally recommended to do increments of 16 wells.
- Fill in the number of cells you will load per transfection (per well)
- The void volume determines the amount of overage to allow for pipetting. The suggested default is 20% which is generally sufficient when using a multichannel pipette and one row of a v-bottom plate as repository for preparing the solution to transfer to the 96-well Nucleocuvette™ Plate.

b) Substrates

Fill in the amount per reaction and stock concentration for the substrate you wish to use. The concentration of pmaxGFP™ Vector control is already entered. For DNA transfections, enter the amount you wish to use per reaction in µg and the concentration of the stock in µg/µl. For siRNA, you may either enter the µg per reaction and stock as µg/µl or the final concentration in a reaction as µM and the stock solution as µM.

GENERAL	Amount	Unit	Default
number of transfections		#	96
cells / transfection		#	
void volume comp. 20µl + X %		%	20

SUBSTRATES		
pmaxGFP™ / reaction		hā
pmaxGFP™ concentration (stock)	0.2	µg/µl
DNA / transfection		hā
DNA concentration (stock)		µg/µl
siRNA / transfection in µg		μg
siRNA concentration (stock)		µg/µl
siRNA / transfection in µM		μΜ
siRNA concentration (stock)		μM

2) Press Calculate to Display the Results in Substrate Volumes, Solutions and Cells

- a) The volume of substrate per well is calculated including 20% overage. This allows you to either prepare individual samples or a master mix in case all samples will get the same substrate.
- b) The solution values will tell the total amount of supplement and Nucleofector™ Solution to mix for all wells and allow for the overage.
- c) The cells value represents the total number of cells to resuspend in the total supplemented Nucleofector™ Solution.

SUBSTRATE VOLUMES pmaxGFP™ (per well) μl pmaxGFP™ (total) μĺ DNA (per well) μl DNA (total) μΙ siRNA (µg-based, per well) μl siRNA (µg-based, total) μl siRNA (µM-based, per well) μΙ siRNA (µM-based, total) μl SOLUTIONS Nucleofector™ Solution μl Supplement μl total volume μΙ CELLS cell number in total volume

3) Rinsing Medium and Cell Culture Post Transfection

These sections can be used to quickly calculate the amount of media to have prewarmed and preplated for post Nucleofection™. Enter the volume to be added per well post Nucleofection™ in the rinsing section and press calculate to get the total prewarmed media needed for adding directly to the Nucleocuvette™ Plate. Enter the Transfer Volume and total plating volume in the Cell culture section to determine the amount to plate prior to beginning Nucleofection™.

RINSING MEDIUM POST TRANSFECT	ION		
medium per well post transfection		μΙ	80
min. vol. of prewarmed medium		ml	
CELL CULTURE POST TRANSFECTION	N		
transfer volume to cell culture		μl	25/50
total cell culture volume		μl	
preplated medium volume per well		μl	

4) Example

transfer volume to cell culture

preplated medium volume per well

total cell culture volume

In the example below, we will be preparing for an experiment with Jurkat cells with 32 samples using siRNA at a final concentration per reaction of $0.5\,\mu\text{M}$ and following the protocol recommendations of 200,000 cells per reaction, 80 μ I media added post Nucleofection, and 50 μ I transferred.

2.0		
32	#	96
200000	#	
20	%	20
	μg	
0.2	µg/µІ	
	μg	
	µg/µl	
	μg	
	µg/µІ	
0.5	μΜ	
20	μΜ	
	ul	
	ш	
No.	μΙ	
	μΙ	
	μΙ	
0.6	μΙ	
19.2	μΙ	
628.2	ul	
	101	
768.0	μl	
-		
76e+6	#	
TION		
80	μl	80
2.56	ml	
	0.2 0.5 0.6 19.2 628.2 139.8 768.0	20 % pg pg/pl pg/pl pg/pl pg/pl pg/pl pg/pl po.5 pM po.5 pM pl

25/50

ul

μΙ

μΙ

50

200

150

Contact Information

North America

Customer Service: 800 638 8174 (toll free)

order.us@lonza.com

Scientific Support: 800 521 0390 (toll free)

scientific.support@lonza.com

Europe

Customer Service: +32 87 321 611

order.europe@lonza.com

Scientific Support: +32 87 321 611 scientific.support.eu@lonza.com

International

Contact your local Lonza distributor
Customer Service: +1 301 898 7025
scientific.support@lonza.com

International Offices

Australia +61 3 9550 0883 Belgium +32 87 321 611 +55 11 2069 8800 Brazil France 0800 91 19 81 (toll free) 0800 182 52 87 (toll free) Germany India +91 40 4123 4000 Ireland 1 800 654 253 (toll free) 800 789 888 (toll free) Italy Japan +81 3 6264 0660 Luxemburg +32 87 321 611 Poland +48 781 120 300 Singapore +65 6521 4379 900 963 298 (toll free) Spain The Netherlands 0800 022 4525 (toll free) United Kingdom 0808 234 97 88 (toll free)

Lonza Cologne GmbH – 50829 Cologne Germany

For research use only. Not for use in diagnostic procedures or for testing or treatment in humans.

Unless otherwise noted, all trademarks herein are marks of the Lonza Group or its affiliates. The information contained herein is believed to be correct and corresponds to the latest state of scientific and technical knowledge. However, no warranty is made, either expressed or implied, regarding its accuracy or the results to be obtained from the use of such information and no warranty is expressed or implied concerning the use of these products. The buyer assumes all risks of use and/or handling. No statement is intended or should be construed as a recommendation to infringe any existing patent.

© Copyright 2012, Lonza Cologne GmbH. All rights reserved. FL-NucCalc 07/12 CD-MN035