

# Protocol for Cryopreservation of Nucleofection™ Competent Cells

Nucleofection™ PLUS Supplements can be used to cryopreserve ready-to-use primary cells or cell lines in Nucleofection™ Solution. For this purpose, the standard supplement must be replaced by Nucleofection™ PLUS Supplement when preparing the respective Nucleofection™ Solution. Once suspended in supplemented Nucleofection™ PLUS Solution cells can be stored in either the liquid nitrogen phase or in the gas phase. We recommend storage in the gas phase in general as liquid nitrogen may seep into the cryovial potentially causing the vial to “burst” upon thawing.

## Freezing Guidelines

### Note:

Cells to be frozen should be free of contamination and in the appropriate phase of growth (see dedicated optimized Nucleofection™ Protocol for further details).

1. Grow the cell line or isolate the primary cell as described in the optimized Nucleofection™ Protocol.
2. **Replace** the standard supplement delivered with the Nucleofection™ Kit by the dedicated Nucleofection™ PLUS Supplement. To find out which Nucleofection™ PLUS Supplement is compatible to a certain Nucleofection™ Kit, please check the list at: [www.lonza.com/n-plus](http://www.lonza.com/n-plus).
3. Allow Nucleofection™ Solution and Nucleofection™ PLUS Supplement to equilibrate at room temperature for 15 minutes.
4. Add the Nucleofection™ PLUS Supplement to the Nucleofection™ Solution.
5. Count an aliquot of the cells and determine cell number.
6. Centrifuge the required number of cells as described in the optimized Nucleofection™ Protocol and carefully discard the supernatant.
7. Re-suspend the cell pellet carefully in appropriate volume of supplemented Nucleofection™ PLUS Solution (20 or 100 µl per planned Nucleofection™ Sample depending on Nucleofection™ Vessel to be used). Accomplish this step at room temperature.
8. Aliquot the cells in cryovials as quick as possible. The cells should not stay in Nucleofection™ PLUS Solution for more than **10 minutes** prior freezing.

### Note:

The size of the cryopreserved cell aliquots should be in accordance with the Nucleofection™ Experiments planned. It is recommended choosing cryovials of a size suited for the total volume of the cryopreserved cell aliquots (recommendation for minimum filling: 25% of the vial volume). Storage of small cell aliquots in huge cryovials bears the risk of water

sublimating to the lid of the cryovial. Due to this phenomenon, the physical parameters of the cell aliquots will change over time which may lead to reduced performance of the cryopreserved cells.

9. Freeze the cell aliquots in a controlled manner by decreasing the temperature by 1°C per minute down to at least to -80°C (recommendation: use a manual device e.g. CoolCell®, Biocision for this purpose). The cells can be stored at -80°C for up to 24 hours prior to the transfer to the cryogenic storage. If you use a controlled rate freezer, decrease the temperature to at least below -125°C before you transfer the vials to the cryogenic storage.
10. Store the Nucleofection™ Competent Cells in the cryogenic storage until they will be used for a Nucleofection™ Experiment.

### Note:

Nucleofection™ Competent Cells are stable for several months to years if stored properly. However, after one year of storage transfection performance may decrease. Therefore it is recommended to run a test transfection before using cells that were stored for a prolonged period of time (more than 12 months).

## Thawing Guidelines

11. Retrieve the cells from the cryogenic storage and transfer them immediately to a 37°C water bath. The cells must thaw quickly (**1 to 2 minutes**) as the temperature range between -50°C and 0°C bears the greatest risk for cellular damage. Thus, passing this temperature range as quickly as possible preserves viability of the competent cells.
12. Gently remove the cells from the vials and transfer them into the appropriate Nucleofection™ Vessels.
13. Mix the cells with substrate of choice by gentle pipetting.
14. Continue with the Nucleofection™ Process as described in the dedicated optimized protocol.

### Note:

Nucleofection™ Competent Cells must be transfected as quickly as possible after thawing and addition of the substrate. The time window mentioned in the cell-type specific Nucleofection™ Protocol does **NOT** apply for cells cryopreserved in Nucleofection™ PLUS Solution. The viability will substantially decrease if thawed cells are kept at room temperature for more than **10 minutes** after thawing.

www.lonza.com  
www.lonza.com/n-plus

## 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: +49 221 99199 400  
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
Brazil	+55 11 5641 3325
France	0800 91 19 81 (toll free)
Germany	0800 182 52 87 (toll free)
India	+91 40 4123 4000
Ireland	1 800 654 253 (toll free)
Italy	800 789 888 (toll free)
Japan	+81 3 6264 0660
Luxemburg	+32 87 321 611
Poland	+48 781 120 300
Singapore	+65 6521 4379
Spain	900 963 298 (toll free)
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.  
The Amaxa™ Nucleofector™ Technology is covered by patent  
and/or patent pending rights owned by Lonza Cologne GmbH.

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.  
P-NucPlusSupp 04/12 CD-KI032

---