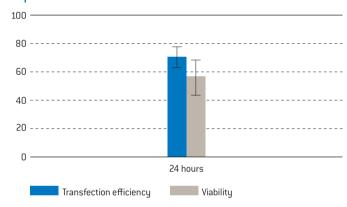


# Amaxa™ HT Nucleofector™ protocol for stimulated human T cells

# Cell description

Stimulated CD3+ human T cells (small, round suspension cells (lymphocyte) are a subpopulation of human peripheral blood mononuclear cells (PBMCs). PBMCs purified from fresh human blood samples treated with anticoagulant or from leucocyte rich buffy coat.

## Example for Nucleofection™ of stimulated human T cells



Transfection efficiency of stimulated human T cells 24 hours post Nucleofection™.  $1\times10^6$  stimulated T cells were transfected with program E0-115-AA (high functionality)using  $0.4~\mu g$  pmaxGFP™ vector. Cells were analyzed 24 hours post Nucleofection™using a FACSCalibur™ with HTS option (Becton Dickinson). Cell viability (CellTiterGlo™Viability Assay, Promega Cat. No.: G7570) is approximately 60 % after 24 hours.

# **Product description**

Recommended kits -P3 Primary Cell HT Nucleofector™ kits

| Cat. No.   | V5SP-3002 |
|--|-----------|
| Size (reactions)                                 | 2×384     |
| P3 primary cell HT Nucleofector™ solution        | 22.5 ml   |
| Supplement                                       | 5 ml      |
| pmaxGFP™ vector (1.0 µg/µl in 10 mM Tris pH 8.0) | 50 μg     |
| 384-well Nucleocuvette <sup>™</sup> plate(s)     | 2         |
| Cat. No.   | V5SP-3010 |

| Cat. No.   | V5SP-3010 |
|--|-----------|
| Size (reactions)                                 | 10×384    |
| P3 primary cell HT Nucleofector™ solution        | 90 ml     |
| Supplement                                       | 20 ml     |
| pmaxGFP™ vector (1.0 μg/μl in 10 mM Tris pH 8.0) | 150 µg    |
| 384-well Nucleocuvette™ plate(s)                 | 10        |

### Storage and stability

Store Nucleofector<sup>™</sup> solution, supplement and pmaxGFP<sup>™</sup> vector at  $4^{\circ}$ C. For long term storage pmaxGFP<sup>™</sup> vector is ideally stored at -20°C. The expiry date is printed on the solution box. Once the Nucleofector<sup>™</sup> supplement is added to the Nucleofector<sup>™</sup> solution it is stable for three months at  $4^{\circ}$ C.

# Note

HT Nucleofector™ solutions can only be used with conductive polymer cuvettes, i.e. in the HT Nucleofector™, the 96-well Shuttle™ device and in the 4D-Nucleofector™ system. They are not compatible with the Nucleofector™ II/2b device.

# Required material

#### Note

Please make sure that the entire supplement is added to the Nucleofector™ solution.

- -HT Nucleofector™
- -Supplemented HT Nucleofector™ solution at room temperature
- -Supplied 384-well Nucleocuvette™ plates
- -Supplied pmaxGFP™ vector, stock solution 1 µg/µl

#### Note

Volume of substrate solution added to each sample should not exceed 10 % of the total reaction volume (2  $\mu$ l for 20  $\mu$ l reactions). For positive control using pmaxGFP<sup>m</sup> vector, please dilute the stock solution to reach the appropriate working concentration.

- Substrate of interest, highly purified, preferably by using endotoxin free Kits; A260 : A280 ratio should be at least 1:8
- 384-well Nucleocuvette™ plates are best pipetted with automated liquid handling systems. If manual pipetting cannot be avoided, please use compatible tips such as: epT.I.P.S.™ (US/CDN: Eppendorf North America, Cat. No. 2491.431, Rest of World: Eppendorf AG, Cat. No. 0030073.266), Matrix TallTips™ (Matrix Technologies Corp., Cat. No. 7281) or LTS Tips (Rainin Instrument, LLC, Cat. No. SR-L10F, SR/SS-L250S, SR/SS-L300S). Please ensure the chosen pipette tips reach the bottom of the 384-well Nucleocuvette™ wells without getting stuck.
- Anti-CD3/anti-CD28 coated HT and 6-well culture plates (see below) or coated culture plates of your choice
- Culture medium: Clonetics™ Lymphocyte Growth Media-3 LGM-3™ for serum-free culture (Lonza, Cat.No. CC-3211) or BioWhittaker™ IMDM media for addition of 10 % serum (Lonza, Cat.No. BE12-722F)
- For isolation: Ficoll-Paque™ Plus (GE Healthcare; Cat. No. 17-1440-03);
   PBS containing 0.5 % (w/v) BSA (PBS/BSA)
- For enrichment (optional): Pan T Cell Isolation Kit II (Miltenyi Biotec;
   Cat. No. 130-091-156) or RosetteSep™ Isolation Kit for human T cells
   (StemCell Technologies, Cat. No 15021)
- For coating of plates (for stimulation): Anti-Human CD3 MAB (OKt 3; eBioscience, Cat. No. 14-0037-82) and Anti-Human CD28 MAB (5E8; Research Diagnostics Inc., Cat. No. 10R-CD28bHUμg/μI); control antibody (purified mlgG(K); BD-Pharmingen, Cat. No. 554 721); antibodies should be diluted in carbonate buffer (32 mM Na2CO3/16 mM NaHCO3) from 100 ng/μI stock solutions directly before use; Immuno™ Plate C96 Maxi Sorp™ (Nunc, Cat. No.: 430 341)
- Prewarm appropriate volume of culture media at 37°C (>95 μl per sample)
- $-\,$  Appropriate number of cells (1×10 $^{\rm 6}$  cells per sample; 5×10 $^{\rm 5}$  cells can be used with slightly reduced transfection efficiency and viability; at even lower cell numbers transfection efficiency and viability are significantly decreased)

# 1. Pre Nucleofection™

#### Note

This protocol is designed for fresh primary human T cells from whole PBMCs. Depending on application T cells can be further enriched (see below).

- Transfection results may be donor-dependent
- For preparation, do not perform protocols using hypo-osmolar buffers.
   This may lead to high cell mortality after Nucleofection™
- For Nucleofection™ of unstimulated T cells, please refer to the optimized protocol for unstimulated human T cells

# Coating of culture plates with anti-CD3 and anti-CD28 antibodies

- 1.1 Incubate each well with 1 ml (for 6-well) or 50 µl (for 96-well; Nunc Immuno™ Plate C96 Maxi Sorp™) of a solution of Anti-Human CD3 MAB at a final concentration of 1 µg/ml and Anti-Human CD28 MAB at a final concentration of 2 µg/ml at 37°C/5 % CO₂ for 5 hours
- 1.2 Wash the wells carefully three times with PBS/BSA

#### **Blood samples**

1.3 Fresh human blood treated with an anticoagulant (e.g. heparin, citrate, ACD-A) or alternatively, leukocyte-enriched buffy coat not older than 8 hours

# **Preparation of PBMC**

- 1.4 Pipet 15 ml Ficoll- Paque™ Plus in a 50 ml conical tube
- 1.5 Overlay Ficoll- Paque™ Plus with 35 ml blood sample and centrifuge at 750×g for 20 minutes at 20°C in a swinging-bucket rotor without brake
- 1.6 Remove the upper layer leaving the mononuclear cell layer undisturbed at the interphase. Carefully transfer the interphase cells (lymphocytes and monocytes) to a new 50 ml conical tube
- 1.7 Add PBS/BSA to 50 ml mark, mix and centrifuge at  $350 \times g$  for 10 minutes at 4°C. Remove the supernatant carefully
- 1.8 Resuspend the cell pellet in 25 ml of PBS/BSA and centrifuge at 160×g for 15 minutes at 4°C. Remove the supernatant carefully
- 1.9 Resuspend the cell pellet in 25 ml PBS/BSA and centrifuge at 300×g for 10 minutes at 4°C. Remove the supernatant carefully
- 1.10 Resuspend cell pellet in 5 ml PBS/BSA and count the cells

#### Note

Purified PBMC may be stored at 4°C overnight in PBS/BSA, but this can cause a significant loss of transfection efficiency.

## Enrichment of T cells (optional)

1.11 Primary human T cells can be further enriched by using Pan T Cell Isolation Kit II (Miltenyi) or RosetteSep™ Isolation Kit for human T cells (StemCell Technologies) according to the manufacturer's protocol

#### Stimulation

1.12 Stimulate the isolated human T cells for 2–3 days prior Nucleofection™ e.g. in plates coated with anti-CD3 antibody and anti-CD28 antibody (please see 1.1–1.2). Seed cells at 5×10<sup>6</sup> cells per ml

# 2. Nucleofection™

## One Nucleofection™ sample contains

- 1×10<sup>6</sup> cells
- 0.4–0.8 μg plasmid DNA (in 1–2 μl  $H_2$ 0 or TE) or 0.4 μg pmaxGFP<sup>™</sup> vector or 30–300 nM siRNA (0.6–6 pmol/sample)
- 20 μl P3 primary cell HT Nucleofector™ solution
- 2.1 Please make sure that the entire supplement is added to the Nucleofector™ solution
- 2.2 Start HT Nucleofector™ Software, verify device connection and upload experimental parameter file (for details please refer to the HT Nucleofector™ manuals)
- 2.3 Select the appropriate HT Nucleofector™ program E0-115-AA
- 2.4 Prepare cell culture plates by filling appropriate number of wells with desired volume of recommended culture media, e.g. 54 µl for one well of a 384-well microtiter plate and pre-incubate/equilibrate plates in a humidified 37°C/5 % CO<sub>2</sub> incubator
- 2.5 Pre-warm an aliquot of culture media to  $37^{\circ}$ C (40  $\mu$ l per sample\* see comments at the end of this chapter)
- 2.6 Prepare 0.4–0.8 µg plasmid DNA or 0.4 µg pmaxGFP™ vector. For siRNA experiments we recommend to start using 30–300 nM siRNA (0.6–6 pmol/sample)
- 2.7 Count the cells and determine cell density
- 2.8 Centrifuge the required number of cells  $(1 \times 10^6)$  cells per sample at 200×g for 10 minutes at room temperature
- 2.9 Resuspend the cell pellet carefully in 20 µl room temperature Nucleofector™ solution per sample

## A: One or several substrates (DNAs or RNAs) in multiples

- Prepare mastermixes by dividing cell suspension according to number of substrates
- Add required amount of substrates to each aliquot (max. 2 µl per sample)

Transfer 20 μl of mastermixes into the wells of the 384-well
 Nucleocuvette™ plate

## B: Multiple substrates (e.g. library transfection)

- Pipette 20  $\mu$ l of cell suspension into each well of a sterile U-bottom 384-well microtiter plate
- Add 2 μl substrates (maximum) to each well

#### Note

It is advisable to pre-dispense each cell suspension into a sterile round bottom 384-well plate or to pipet from a pipetting reservoir for multichannel pipettes. Use a liquid handling system or manual pipette with suitable pipette tips. As leaving cells in HT Nucleofector™ solution for extended periods of time may lead to reduced transfection efficiency and viability it is important to work as quickly as possible. Avoid air bubbles while pipetting.

- 2.10 Briefly shake 384-well Nucleocuvette™ plate with an appropriate microtiter plate shaker to make sure the sample covers the bottom and sides of the well with no air bubbles. Alternatively, thoroughly tap the 384-well Nucleocuvette™ plate
- 2.11 Place 384-well Nucleocuvette™ plate with closed lid onto the carousel of the Plate Handler. Well "A1" must be in upper left position
- 2.12 Start HT Nucleofection™ process by clicking "Start" in the HT Nucleofector™ Software
- 2.13 After run completion, carefully remove the 384-well Nucleocuvette™ plate from the carousel
- 2.14 Resuspend cells with desired volume of pre-warmed media (maximum cuvette volume 60 µl). Mix cells by gently pipetting up and down two to three times. Recommendation for 384-well Nucleocuvette™ plates: Resuspend cells in 40 µl of pre-warmed media\*
- 2.15 Plate desired amount of cells in culture system of your choice. Recommendation for 384-well cell culture plates: Transfer 6  $\mu$ l of resuspended cells to 54  $\mu$ l pre-warmed media prepared in 384-well cell culture plates\*

#### \* Note

The indicated cell numbers and volumes have been found to produce optimal Nucleofection™ results in most cases, however, depending on your specific needs you may wish to test an extended range of cell numbers. cell numbers and volumes can be adapted such that fewer cells are transferred or duplicate plates can be seeded.

# 3. Post Nucleofection™

- 3.1 Incubate the cells in a humidified 37°C/5 % CO<sub>2</sub> incubator until analysis. Gene expression or down regulation, respectively, is often detectable after only 4–8 hours. If this is not the case, the incubation period may be prolonged to 24 or 48 hours
- 3.2 Culture stimulated T cells post Nucleofection™ in plates coated with anti-CD3 antibody and anti-CD28 antibody (see chapter 1)

# Additional information

# Up-to-date List of all Nucleofector™ references

www.lonza.com/nucleofection-citations

# Technical assistance and scientific support

#### USA/Canada

Tel 800 521 0390 (toll-free) Fax 301 845 8338 scientific.support@lonza.com

Europe and Rest of World

Tel +49 221 99199 400 Fax +49 221 99199 499 scientific.support.eu@lonza.com

#### www.lonza.com

#### Lonza Cologne GmbH-50829 Cologne, Germany

Please note that the Amaxa" Nucleofector" technology is not intended to be used for diagnostic purposes or for testing or treatment in humans.

The Nucleofector" technology, comprising Nucleofection" process, Nucleofector" device, 4D Nucleofector" device, HT Nucleofector device, Nucleofector device, Nucleofector solutions and 96-well Nucleocuvette plates and modules is covered by patent and/or patent-pending rights owned by Lonza Cologne GmbH.

Amaxa, Nucleofector, Nucleofection, maxGFP, Reagent Pack, HCM and BulletKit are either registered trademarks or trademarks of the Lonza Group or its affiliates.

TallTips are a registered trademark of Matrix Technologies Corporation.

Other product and company names mentioned herein are the trademarks of their respective owners.

This kit contains a proprietary nucleic acid coding for a proprietary copepod fluorescent protein intended to be used as a positive control with this Lonza product only. Any use of the proprietary nucleic acid or protein other than as a positive control with this Lonza product is strictly prohibited. USE IN ANY OTHER APPLICATION REQUIRES A LICENSE FROM EVROGEN. To obtain such a license, please contact Evrogen at license@evrogen.com.

The CMV promoter is covered under U.S. Patents 5,168,062 and 5,385,839 and its use is permitted for research purposes only. Any other use of the CMV promoter requires a license from the University of lowa Research Foundation, 214 Technology Innovation Center, Iowa City, IA 52242.

The use of this product in conjunction with materials or methods of third parties may require a license by a third party.

User shall be fully responsible for determining whether and from which third party it requires such license and for the obtainment of such license.

No statement is intended or should be construed as a recommendation to infringe any existing patent.

© Copyright 2011, Lonza Cologne GmbH. All rights reserved DSSP-3007\_2011-10