

## Human CD4<sup>+</sup>CD25<sup>+</sup> regulatory T cell isolation, *in vitro* expansion and analysis

### *In vitro* human regulatory T cell expansion

#### Introduction

Regulatory T (Treg) cells are a subpopulation of T cells. Basic as well as clinical research has focused on the study of Treg cells for years due to their unique function of suppressing immune responses. As Treg cells represent a small cell population, many downstream applications require an expansion of Treg cells prior to analysis. However, Treg cells have been described as hypoproliferative in response to polyclonal stimulation and interleukin 2 (IL-2) *in vitro*. Thus, *in vitro* cultivation of Treg cells often results in low expansion rates or low frequencies of FoxP3<sup>+</sup> Treg cells due to overgrowth of conventional T cells.

To overcome this problem, we have developed the Treg Expansion Kit, human, which is designed to efficiently expand Treg cells that maintain FoxP3 expression after isolation with the CD4<sup>+</sup>CD25<sup>+</sup> Regulatory T Cell Isolation Kit, human.

This application note includes a complete protocol for your Treg cell workflow, describing human Treg cell isolation from peripheral blood mononuclear cells (PBMCs), their *in vitro* expansion and subsequent flow cytometry analysis.

#### Workflow

<b>Day 0</b>	Treg cell isolation  Flow cytometry analysis
<b>Day 1</b>	Addition of 100 µL expansion medium
<b>Day 3-5</b>	Split cells or replace 100 µL medium with fresh expansion medium
<b>Day 5-8</b>	Transfer Treg cells into 24-well plate or split cells and use a new 96-well plate
<b>Day 14</b>	Stop Treg expansion and analyze Treg cells by flow cytometry OR Restimulate Treg cells
<b>Day 21</b>	Stop Treg cell expansion  Flow cytometry analysis

## Materials

### Treg cell isolation from PBMCs

#### Treg cell isolation kit

- CD4<sup>+</sup>CD25<sup>+</sup> Regulatory T Cell Isolation Kit, human (# 130-091-301)

#### Separation buffer

- 1x PBS (pH 7.2), BSA, EDTA
- Alternative: 6x 1.5 L autoMACS<sup>®</sup> Running Buffer – MACS<sup>®</sup> Separation Buffer (# 130-091-221)

#### Magnetic separation

- LD Columns, 1 for 10<sup>8</sup> labeled cells (# 130-042-901)
- MS Columns, 2x per isolation (# 130-042-201)
- MidiMACS<sup>™</sup> Separator (Separator for LD Column), (# 130-042-302)
- MiniMACS<sup>™</sup> Separator (Separator for MS Column), (# 130-042-102)
- MACS<sup>®</sup> MultiStand (# 130-042-303)
- (Optional) Pre-Separation Filters, 30 μm (# 130-041-407)

### Treg cell *in vitro* expansion

#### Expansion medium

- TexMACS<sup>™</sup> Medium, research grade (500 mL), (# 130-097-196)
- Human IL-2 IS, premium grade (50 μg) (# 130-097-745)
- human AB Serum
- β-Mercaptoethanol (β-ME)
- Penicillin/Streptomycin (PenStrep)
- (Optional) Rapamycin

#### Expansion process

- Treg Expansion Kit, human: 2 mL for stimulation of 1x10<sup>7</sup> Treg cells (# 130-095-345); 2x2 mL for stimulation of 2x10<sup>7</sup> Treg cells (# 130-095-353)
- 96-well round-bottom plate
- (Optional) 24-well plate

#### Bead removal (optional)

- MACSiMAG<sup>™</sup> Separator (# 130-092-168)

### Staining of Treg cells for flow cytometry analysis

- Treg Detection Kit (CD4/CD25/FoxP3) (PE), human (# 130-094-163)

OR

- Treg Detection Kit (CD4/CD25/FoxP3) (APC), human (# 130-094-158)
- (Optional): CD127- PE-Vio 770 (# 130-109-516) or CD127- PE-Vio 615, human (# 130-107-513)

## Material preparation

### Treg cell isolation from PBMCs

#### Separation buffer

- 1x PBS (pH 7.2)  
+ 0.5% BSA  
+ 2 mM EDTA

### Treg cell *in vitro* expansion

#### Expansion medium

The expansion medium can be stored for up to 10 days under sterile conditions at 4°C.

- 500 mL TexMACS<sup>™</sup> Medium  
+ 500 IU/mL Human IL-2 IS, premium grade  
+ 5% human AB Serum  
+ 0.01 mM β-ME  
+ 1% PenStrep  
+ (Optional) 100 nmol/L Rapamycin

▲ **Note:** Rapamycin has been shown to increase the frequency of FoxP3<sup>+</sup> Treg cell during expansion (Battaglia, M. *et al.* (2005) *Rapamycin selectively expands CD4+CD25+FoxP3+ regulatory T cells. Blood 105: 4743–4748*).

### Flow cytometry analysis of Treg cell expansion (intracellular staining)

#### FoxP3 staining buffer

(components provided in Treg Detection Kit)

▲ Always prepare reagents freshly on the same day that intracellular FoxP3 staining is performed.

▲ The required total buffer volumes should be calculated beforehand; volumes will depend on the number of cells to be analyzed as well as the number of tests to be performed.

### Example:

For 1 sample with  $1 \times 10^6$  Treg cells: prepare 1 mL of fixation/permeabilization solution and 4 mL of 1× Permeabilization Buffer

#### I) Fixation/permeabilization solution

- Dilute Fixation/Permeabilization Solution 1 (e.g. 0.25 mL) 1:4 with Fixation/Permeabilization Solution 2 (e.g. 0.75 mL).

#### II) 1× Permeabilization Buffer

- Dilute the 10× Permeabilization Buffer (e.g. 1 mL) 1:10 with deionized or distilled water (e.g. 9 mL).

## Methods

### 1. Generation of PBMCs

When working with anticoagulated peripheral blood or buffy coat, PBMCs should be isolated by density gradient centrifugation, for example, using Ficoll-Paque™ according to the manufacturer's instructions. Collect a sample of PBMCs for subsequent flow cytometry analysis.

### 2. Isolation of CD4<sup>+</sup>CD25<sup>+</sup> Treg cells using the CD4<sup>+</sup>CD25<sup>+</sup> Regulatory T Cell Isolation Kit, human

#### 2.1 Magnetic labeling of non-CD4<sup>+</sup> cells

▲ Volumes for magnetic labeling given below are for up to  $10^7$  total cells. When working with fewer than  $10^7$  cells, use the same volumes as indicated. When working with higher cell numbers, scale up all reagent volumes and total volumes accordingly (e.g. for  $2 \times 10^7$  total cells, use twice the volume of all indicated reagent volumes and total volumes)

1. Determine cell number.
2. Centrifuge cell suspension at 300×g for 10 minutes. Discard supernatant.
3. Resuspend cell pellet in 90 μL of separation buffer per  $10^7$  total cells.
4. Add 10 μL of CD4<sup>+</sup> T Cell Biotin-Antibody Cocktail per  $10^7$  total cells.

5. Mix well and incubate for 5 minutes in the refrigerator (2–8 °C).

6. Add 20 μL of Anti-Biotin MicroBeads per  $10^7$  total cells.

7. Mix well and incubate for 10 minutes in the refrigerator (2–8 °C).

8. Adjust the volume to a minimum of 500 μL of separation buffer.

▲ **Note:** Resuspend up to  $10^8$  cells in 500 μL of separation buffer. For higher cell numbers, scale up buffer volume accordingly.

#### 2.2 Magnetic depletion of non-CD4<sup>+</sup> cells

The first step of Treg cell isolation is a depletion of non-CD4<sup>+</sup> cells. Here, an LD Column is used, which has a capacity of  $10^8$  labeled cells and  $5 \times 10^8$  total cells.

▲ **Note:** When using the CD4<sup>+</sup>CD25<sup>+</sup> Regulatory T cell Isolation Kit, human, we do not recommend to process more than  $1.3 \times 10^8$  total cells on an LD Column. When exceeding this cell number, it is strongly recommended to split the sample and use additional LD Columns.

1. Place LD Column(s) in the magnetic field of a MidiMACS™ Separator.

2. Prepare column by rinsing with 2 mL of separation buffer. Always wait until the column reservoir is empty before proceeding to the next step.

3. Apply cell suspension onto the column.

4. Collect unlabeled cells that pass through and wash column twice with 1 mL of separation buffer each. **Collect total effluent; this is the unlabeled pre-enriched CD4<sup>+</sup> cell fraction which is needed for further Treg cell isolation.**

5. Determine cell number.

#### 2.3. Magnetic labeling of CD25<sup>+</sup> cells

▲ Volumes for magnetic labeling given below are for an initial starting cell number of up to  $10^7$  total cells. For higher initial cell numbers, scale up all volumes accordingly.

1. Centrifuge cell suspension at 300×g for 10 minutes. Discard supernatant.

2. Resuspend cell pellet in 90  $\mu\text{L}$  of separation buffer per  $10^7$  total cells.

3. Add 10  $\mu\text{L}$  of CD25 MicroBeads II per  $10^7$  total cells.

4. Mix well and incubate for 15 minutes in the refrigerator (2–8  $^{\circ}\text{C}$ ).

5. Wash cells by adding 1–2 mL of separation buffer and centrifuge at  $300\times g$  for 10 minutes. Discard supernatant.

6. Resuspend up to  $10^8$  cells in 500  $\mu\text{L}$  of separation buffer.

▲ **Note:** For higher cell numbers, scale up buffer volume accordingly.

#### 2.4. Magnetic separation of CD25<sup>+</sup> cells

The second step during Treg cell isolation is a positive selection of CD25<sup>+</sup> cells. Here, two consecutive MS Columns are used, with a capacity of  $10^7$  labeled cells. To not exceed the capacity, it is recommended to determine the frequency of CD25<sup>+</sup> cells in your cell suspension by flow cytometry beforehand.

1. Place an MS Column in the magnetic field of a MiniMACS™ Separator.

2. Prepare column by rinsing with 500  $\mu\text{L}$  of separation buffer.

3. Apply cell suspension onto the column. Collect flow-through containing unlabeled cells.

4. Wash column 3 times with 500  $\mu\text{L}$  of separation buffer each. Collect unlabeled cells that pass through and combine with the effluent from step 3.

5. Remove column from the separator and place it on a suitable collection tube.

6. Pipette 1 mL of separation buffer onto the column. Immediately flush out magnetically labeled cells by firmly pushing the plunger into the column.

7. To increase purity of CD4<sup>+</sup>CD25<sup>+</sup> cells, the eluted fraction can be enriched over a second MS Column (recommended). Repeat the

magnetic separation procedure described in steps 1 to 6 using a new MS Column. **The isolated Treg cells are now ready-to-use for *in vitro* expansion.**

To assess the purity and to determine the frequency of FoxP3-expressing cells as start population for expansion, a surface and intracellular flow cytometry staining of the isolated Treg cells must be performed. The purity is determined by surface staining of CD4 and CD25. Furthermore, for the final analysis of Treg expansion, it is also important to determine:

- a) The initial cell number
- b) The initial cell number and frequency of FoxP3-expressing CD4<sup>+</sup> cells

#### 3. Surface and intracellular flow cytometry staining of Treg cells using the Treg Detection Kit

▲ Volumes given below are **for up to  $10^6$**  nucleated cells. When working with fewer than  $10^6$  cells, use the same volumes as indicated **except for the fixation step** due to the impact on cell morphology. When working with higher cell numbers, scale up all reagent volumes and total volumes, accordingly (e.g. for  $2\times 10^6$  nucleated cells, use twice the volume of all indicated reagent volumes and total volumes).

##### 3.1 Surface staining of Treg cells with CD4 and CD25 antibodies

1. Determine cell number.

2. Centrifuge cell suspension at  $300\times g$  for 10 minutes. Discard supernatant.

3. Resuspend up to  $10^6$  nucleated cells per 90  $\mu\text{L}$  of separation buffer.

4. Add 10  $\mu\text{L}$  of CD4-FITC and 20  $\mu\text{L}$  of CD25 antibody (either PE or APC conjugated).

5. Mix well and incubate for 10 minutes in the dark in the refrigerator (2–8  $^{\circ}\text{C}$ ).

6. Wash cells by adding 1–2 mL of separation buffer per  $10^6$  cells and centrifuge at  $300\times g$  for 5 minutes at 4  $^{\circ}\text{C}$ . Discard supernatant

7. Proceed immediately to 3.2.

### 3.2 Intracellular staining with the Anti-FoxP3 antibody

1. Resuspend  $10^6$  nucleated cells in 1 mL cold, freshly prepared fixation/permeabilization solution (for details refer to material preparation section). When working with fewer cells, scale down the buffer volume accordingly (e.g. for  $5 \times 10^5$  nucleated cells use 500  $\mu$ L of fixation/permeabilization solution).

2. Mix well and incubate for 30 minutes in the dark in the refrigerator (2–8 °C).

3. Wash cells by adding 1–2 mL of cold separation buffer per  $10^6$  cells and centrifuge at  $300 \times g$  for 5 minutes at 4 °C. Discard supernatant.

4. Wash cells by adding 1–2 mL of cold 1 $\times$  Permeabilization Buffer (for details refer to material preparation section) per  $10^6$  cells and centrifuge at  $300 \times g$  for 5 minutes at 4 °C. Discard supernatant.

5. Resuspend up to  $10^6$  nucleated cells in 80  $\mu$ L of cold 1 $\times$  Permeabilization Buffer.

6. Add 20  $\mu$ L of FcR Blocking Reagent.

7. Mix well and incubate for 5 minutes in the refrigerator (2–8 °C).

8. Add 10  $\mu$ L of the Anti-FoxP3 antibody (either APC or PE conjugated).

9. Mix well and incubate for 30 minutes in the dark in the refrigerator (2–8 °C).

10. Wash cells by adding 1–2 mL of cold 1 $\times$  Permeabilization Buffer per  $10^6$  cells and centrifuge at  $300 \times g$  for 5 minutes at 4 °C. Discard supernatant.

11. Resuspend cell pellet in a suitable amount of separation buffer for analysis by flow cytometry.

▲ **Note:** Due to fixation and permeabilization, cells can be smaller than viable cells. Thus, FSC/SSC settings of the flow cytometer might need to be adjusted.

### 3.3. Flow cytometry analysis

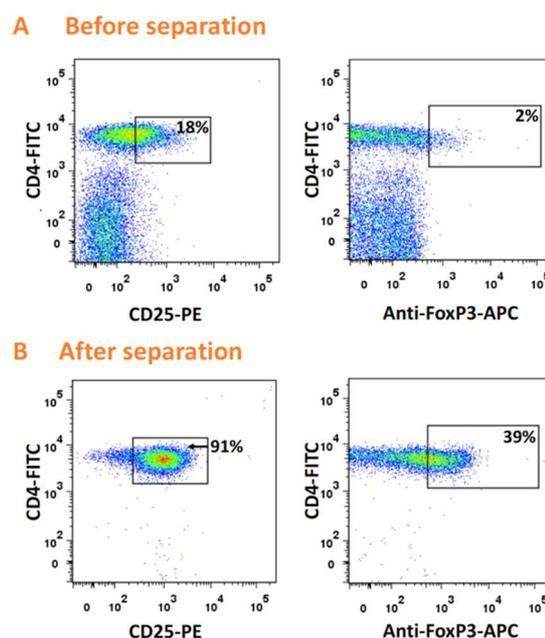
To assess the purity of the isolated Treg cells and the initial frequency of FoxP3<sup>+</sup> CD4<sup>+</sup> cells, the cells were analyzed by flow cytometry. The analysis should be performed with the cell sample taken in section 1 (before separation) and cell sample after isolation (after separation).

1. Lymphocytes were identified by FSC/SSC gating (data not shown).

2. Dead cells were excluded from the analysis by using the live/dead cell exclusion marker PI (data not shown).

3. Lymphocytes were further analyzed for their expression of CD4 (y-axis) and CD25 (x-axis) to assess the purity of CD4<sup>+</sup>CD25<sup>+</sup> Treg cells (fig. 1 A and B, left plots).

4. Lymphocytes were further analyzed for their expression of CD4 (y-axis) and FoxP3 (x-axis) to assess the initial frequency of FoxP3-expressing CD4<sup>+</sup> cells among lymphocytes (fig. 1 A and B, right plots).



**Figure 1: Flow cytometry analysis of Treg cells.**

CD4<sup>+</sup>CD25<sup>+</sup>regulatory T cells were isolated from human PBMCs by using the CD4<sup>+</sup>CD25<sup>+</sup> Regulatory T Cell Isolation Kit, human. The cells were fluorescently stained with CD4-FITC, CD25-PE and Anti-FoxP3-APC before (A) and after (B) separation.

#### 4. *In vitro* expansion of regulatory T cells using the Treg Expansion Kit

Prepare the expansion medium as described in section “Material preparation”. The Treg Expansion Kit is based on MACSiBead™ Particles preloaded with CD3 and CD28 antibodies. Best expansion of Treg cells is accomplished by using MACSiBead Particles and Treg cells at a bead-to-cell ratio of 4:1.

##### 4.1. Preparation of Treg cells

1. Determine the total number of Treg cells.
2. Wash Treg cells by adding 5–10 volumes expansion medium to the cells and centrifuge at 300×g for 10 minutes. Discard supernatant.
3. Resuspend cells at a concentration of  $1 \times 10^6$  cells/mL in expansion medium. Pipet 100  $\mu$ L ( $1 \times 10^5$  Treg cells) in each well of a 96-well round-bottom plate.

##### 4.2. Preparation of CD3/CD28 MACSiBead™ Particles

▲ The CD3/CD28 MACSiBead™ Particles have a concentration of  $2 \times 10^7$  beads/mL and should be used at a bead-to-cell ratio of 4:1.

▲ MACSiBead Particles are bigger in size than MACS® MicroBeads and sediment rapidly. It is therefore mandatory to bring the MACSiBead Particles into suspension by vortexing prior to use.

1. Resuspend CD3/CD28 MACSiBead Particles thoroughly by vortexing and transfer 4 times the amount of Treg cells to a suitable tube. (e.g. for expansion of  $1 \times 10^6$  Treg cells, transfer  $4 \times 10^6$  (=200  $\mu$ L) of MACSiBead Particles).
2. Wash MACSiBead Particles by adding 300–600  $\mu$ L of culture medium and centrifuge at 300×g for 5 minutes. Aspirate supernatant completely.
3. Resuspend CD3/CD28 MACSiBead Particles in expansion medium to get a final concentration of  $2 \times 10^7$  beads/mL (e.g. when

$4 \times 10^6$  of MACSiBead Particles are used, resuspend in 200  $\mu$ L of expansion medium).

##### 4.3. Co-cultivation of Treg cells and MACSiBead Particles

1. Add 20  $\mu$ L of CD3/CD28 MACSiBead Particles to each well containing the Treg cells. As the MACSiBead Particles sediment rapidly, the bead suspension should be vortexed from time to time prior addition to the well.
2. The final volume of each well is now 120  $\mu$ L. Let the co-culture of Treg cells and MACSiBead Particles rest for one day at 37°C and 5–7% CO<sub>2</sub>.

##### 4.4. Expansion process

**Day 1:** Add 100  $\mu$ L of expansion medium to each well.

**Day 3-5** (depending on medium color): Either split Treg cells (including MACSiBead Particles) at a ratio of 1:2 and resuspend cells in 200  $\mu$ L of fresh expansion medium or carefully aspirate 100  $\mu$ L of old medium and add 100  $\mu$ L of fresh expansion medium.

**Day 5-8:** Either transfer Treg cells (including MACSiBead Particles) to a 24-well plate in higher volumes, e.g.,  $5 \times 10^5$  Treg cells in 500  $\mu$ L, or split Treg cells (including MACSiBead Particles) at a ratio of 1:2 and resuspend cells in 200  $\mu$ L of fresh expansion medium.

**Day 14:** The Treg expansion process can be stopped here or proceeded by restimulation of Treg cells. If you want to stop Treg expansion here, determine the total cell number and the frequency of CD4<sup>+</sup>FoxP3-expressing cells by performing a surface and intracellular flow cytometry staining to assess the result of expansion. For details refer to section 3. If you want to proceed with restimulation the CD3/CD28 MACSiBead Particles may be removed prior restimulation. For details refer to section 5.

##### 5. Bead removal and restimulation

Removal of MACSiBead Particles may be required before restimulation with MACSiBead Particles. In addition, bead removal is required before magnetic separation of cells with MACS

MicroBeads or before stimulation with different agents or antigens.

### 5.1. Removal of MACSiBead Particles

1. Harvest cells, transfer to an appropriate tube and wash once with separation buffer.

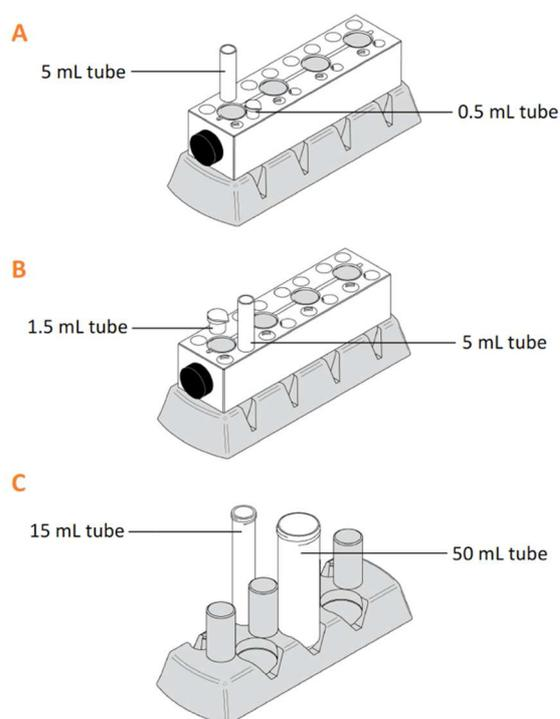
2. Determine cell number and resuspend cells in separation buffer at a density of up to  $2 \times 10^7$  cells/mL and vortex thoroughly.

3. Place the tube in the magnetic field of the MACSiMAG™ Separator (fig. 2). Detailed instructions on how to use the MACSiMAG Separator can be found in the MACSiMAG Separator datasheet.

4. Allow the MACSiBead Particles to adhere to the wall of the tube:

- 5 mL tubes: 2 minutes
- 15 mL or 50 mL tubes: 4 minutes

5. Retaining the tube in the magnet, carefully remove the supernatant containing the MACSiBead Particles-depleted cells and place in a new tube.



**Figure 2: The MACSiMAG Separator.** (A) MACSiMAG Separator with tube rack positioned for tubes of 5 mL or 0.5 mL in size. (B) MACSiMAG Separator with tube rack positioned for tubes of 1.5 mL or 5 mL in size. (C) MACSiMAG Separator with 15 mL and 50 mL tube.

6. Remove the tube from the separator and add separation buffer to reach the same volume as before.

7. Vortex sample, replace tube in the MACSiMAG Separator and repeat steps 4–5.

8. Collected cells can now be further processed as required, e.g., for restimulation with MACSiBead Particles.

### 5.2. Restimulation of Treg cells

▲ Restimulation is performed with a bead-to-cell ratio of 4:1.

Depending on Treg number and resulting volumes, the restimulation can either be performed in a 96-well plate or in a 24-well plate. If you prefer to restimulate Treg cells in a 96-well-plate, refer to section 4 and follow the protocol. Alternatively, restimulation may also be performed in 24-well plates in higher volumes, for example,  $5 \times 10^5$  Treg cells in 500  $\mu$ L of expansion medium and 100  $\mu$ L of CD3/CD28 MACSiBead Particles (=  $2 \times 10^6$  particles).

**Day 15:** Add 100  $\mu$ L of expansion medium per well when restimulation is performed in a 96-well plate, or 500  $\mu$ L of expansion medium per well when restimulation is performed in a 24-well plate.

**Day 17-19:** (depending on medium color): Either split Treg cells again (including the MACSiBead Particles) at a ratio of 1:2 and fill the well to 1 mL with fresh expansion medium or carefully aspirate 500  $\mu$ L old medium and add 500  $\mu$ L fresh expansion medium.

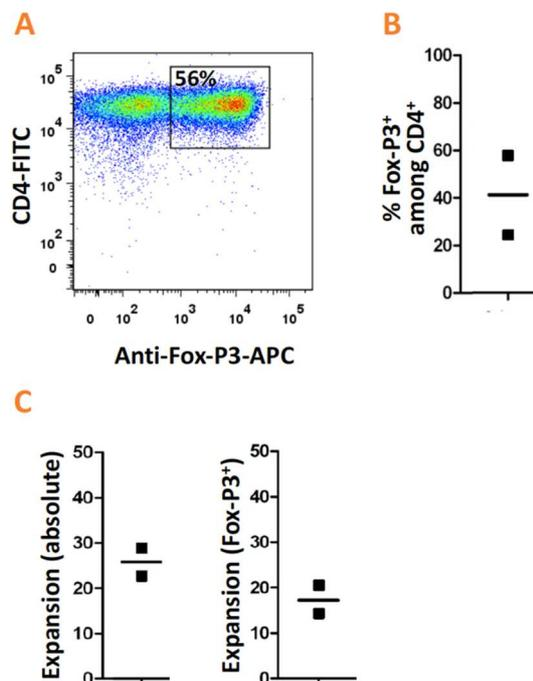
**Day 21:** The Treg expansion process should be stopped after 21 days. To assess the result of expansion, determine the total cell number and the frequency of FoxP3-expressing cells, by performing a surface and intracellular flow cytometry staining. For details see section 3.

▲ **Note:** Bead removal is required before certain downstream applications, i.e. magnetic separation of cells with MACS MicroBeads or before stimulation with different agents or antigens. For details refer to section 5.1.

### 5.3. Flow cytometry analysis

To determine the frequency of FoxP3-positive CD4<sup>+</sup> cells after expansion, the cells were analyzed by flow cytometry.

1. Lymphocytes were identified according to FSC and SSC (data not shown).
2. Dead cells were excluded from the analysis by using the live/dead cell exclusion marker PI (data not shown).
3. Lymphocytes were further analyzed for their expression of CD4 (y-axis) and FoxP3 (x-axis) to assess the frequency of FoxP3-expressing CD4<sup>+</sup> cells among lymphocytes (fig. 3A). Please note that CD4 is used here to identify Treg cells as CD25 is not an appropriate Treg marker after expansion as it might also be upregulated on activated T cells.



**Figure 3: Analysis of Treg expansion.** CD4<sup>+</sup>CD25<sup>+</sup> regulatory T cells were isolated from human PBMCs by using the CD4<sup>+</sup>CD25<sup>+</sup> Regulatory T Cell Isolation Kit, human and expanded for 21 days with the Treg Expansion Kit. Afterwards, the cells were fluorescently stained with CD4-FITC and FoxP3-APC antibodies. (A) Representative FACS plot of expanded Treg cells and (B) frequency of FoxP3<sup>+</sup> Treg cells among CD4<sup>+</sup> cells after expansion from two experiments. (C) Fold-expansion of total cells (left) and FoxP3<sup>+</sup> cells (right) (Fold-expansion comparing the initial and the final cell number). Each dot represents one experiment.

► [miltenyibiotec.com](http://miltenyibiotec.com)

Miltenyi Biotec provides products and services worldwide. Visit [www.miltenyibiotec.com/local](http://www.miltenyibiotec.com/local) to find your nearest Miltenyi Biotec contact.

Unless otherwise specifically indicated, Miltenyi Biotec products and services are for research use only and not for therapeutic or diagnostic use. MACS, the MACS logo, MidiMACS, MiniMACS, TexMACS, MACSiBead, and VioBlue are registered trademarks or trademarks of Miltenyi Biotec GmbH and/or its affiliates in various countries worldwide. All other trademarks mentioned in this document are the property of their respective owners and are used for identification purposes only. Copyright © 2017 Miltenyi Biotec GmbH and/or its affiliates. All rights reserved.