

September 2012

## Simultaneous Measurement of Mouse Foxp3 and Ki-67 in Cultured Splenocytes

### Materials and Reagents

| Full Name  | Short Name              | Catalog Number |
|--|-------------------------|----------------|
| BD Falcon™ polystyrene 12 x 75-mm tubes                    | sample tubes            | 352058         |
| BD Falcon™ 75-cm <sup>2</sup> cell culture flask           | culture flask           | 353136         |
| BD Falcon™ conical tubes with screw cap                    | conical tube            | 358206         |
| RPMI-1640 containing 10% FCS                               | medium                  |                |
| Sterile 1X PBS   | 1X PBS                  |                |
| BD Pharmingen™ Mouse Foxp3 buffer set                      | Foxp3 Buffer Set        | 560409         |
| BD Pharm Lyse™ solution                                    | Lysing Solution         | 555899         |
| BD Pharmingen™ Stain Buffer (BSA)                          | Stain Buffer            | 554657         |
| BD Cytotfix/Cytoperm™ Fixation/Permeabilization Solution   | Fix/Perm Solution       | 554722         |
| BD Cytotfix/Cytoperm™ Perm/Wash™ Buffer                    | Perm/Wash Buffer        | 554723         |
| NA/LE anti-mouse CD3e, clone 145-2C11                      | Anti-CD3                | 553057         |
| NA/LE anti-mouse CD28, clone 37.51                         | Anti-CD28               | 553294         |
| Alexa Fluor® 488 Rat Anti-mouse CD4, clone RM4-5           | CD4- Alexa Fluor® 488   | 557667         |
| PerCP-Cy™5.5 Rat Anti-Mouse CD25, clone PC61               | CD25-PerCP-Cy5.5        | 551071         |
| Alexa Fluor® 647 Rat anti-Mouse FoxPp3, clone MF23         | Foxp3- Alexa Fluor® 647 | 560401         |
| PE Mouse Anti-Human Ki-67 Set                              | Ki-67-PE                | 556027         |
| Anti-Rat Ig, κ/Negative Control Compensation Particles Set | Compensation beads      | 552844         |

### Procedural Notes

- This assay enables the simultaneous measurement of cellular proliferation (using the Ki-67 marker) and cell surface markers in Foxp3-positive cells.
- In this protocol, the cells were prepared from mouse spleen Balb/c mice (female). The mice were sacrificed, spleen was harvested, and the red blood cells were lysed with Lysing Solution. The splenocytes were added to a culture flask at  $1 \times 10^6$  cells/mL of medium.
- Do not store diluted Fixation Buffer and Permeabilization Buffer prepared from the Foxp3 Buffer Set. Prepare them fresh every time.
- Determine the appropriate amount of the Rat Anti-mouse Foxp3 antibody by titration.

### Procedure: Stimulation

1. Coat a sterile culture flask with anti-CD3 at a concentration of 10 µg/mL in sterile 1X PBS. Incubate overnight at room temperature (RT).
2. The next day, discard the anti-CD3 and wash once with 2 mL of medium.
3. Prepare splenocytes from mouse spleen. Lyse the red blood cells for 2 minutes at RT, using 2 mL of Lysing Solution for each spleen.
4. Add 18 mL of RPMI and centrifuge at 250g for 5 to 10 minutes. Remove the supernatant.
5. Resuspend the cells and plate at  $1 \times 10^6$  cells/mL of medium onto a culture flask.
6. Add anti-CD28 at a concentration of 2 µg/mL and incubate for 72 hours at 37°C in 5% CO<sub>2</sub>.
7. In parallel, prepare a mock treated flask using the same protocol but excluding the stimulation antibodies.
8. Harvest the cells after 72 hours. Wash the cells two times with Stain Buffer.
9. Pellet the cells by centrifugation at 250g for 10 minutes. Remove the supernatant.
10. Mix the pellet to loosen the cells. Resuspend the cells at  $10 \times 10^6$  cells/mL of Stain Buffer.
11. Proceed to the Cell Preparation and Foxp3 Staining procedure.



September 2012

## Simultaneous Measurement of Mouse Foxp3 and Ki-67 in Cultured Splenocytes

### Procedure: Buffer Preparation (from Foxp3 Buffer Set)

- Prepare 1X Fixation Buffer.
  1. Dilute 1 part Fixation Concentrate with 19 parts 1X PBS.
  2. Prepare 2 mL for each test ( $1 \times 10^6$  cells).
  3. Place on ice.
- Prepare 1X Permeabilization Buffer.
  1. Dilute 1 part Permeabilization Concentrate with 4 parts 1X PBS.
  2. Prepare 4 mL for each test ( $1 \times 10^6$  cells).
  3. Pre-warm to 37°C before use.

### Procedure: Cell Preparation and Foxp3 Staining

1. Aliquot the appropriate amount of selected surface staining reagents to sample tubes.
2. Add 100  $\mu$ L of cells to each sample tube and mix well. Incubate for 20 minutes at RT in the dark.
3. Wash the cells with 2 mL of Stain Buffer. Pellet the cells by centrifugation at 250g for 10 minutes.
4. Remove the supernatant. Mix the pellet to loosen the cells.
5. Fix the cells with 2 mL of freshly prepared cold 1X Fixation Buffer. Mix well. Incubate for 30 minutes at 4°C in the dark.
6. Pellet the cells by centrifugation at 500g for 5 minutes. Remove the Fixation Buffer.
7. Wash the cells with 2 mL of prepared pre-warmed 1X Permeabilization Buffer.
8. Pellet the cells by centrifugation at 500g for 5 minutes.
9. Remove the supernatant. Mix the pellet to loosen the cells.
10. Permeabilize the cells by adding 2 mL of freshly prepared pre-warmed 1X Permeabilization Buffer. Incubate for 30 minutes at 37°C in the dark.
11. Pellet the cells by centrifugation at 500g for 5 minutes. Remove the Permeabilization Buffer.
12. Wash the cells with 2 mL of Stain Buffer. Pellet the cells by centrifugation at 500g for 10 minutes. Remove the supernatant, leaving about 100  $\mu$ L in the tube. Mix the pellet to loosen the cells.
13. Add 20  $\mu$ L of Rat Anti-Mouse Foxp3 antibody (dilute the appropriate amount in Stain Buffer) to each sample tube and mix well. Incubate for 20 minutes at RT in the dark.
14. Wash the cells twice with 2 mL of Stain Buffer. Pellet the cells by centrifugation at 500g for 10 minutes. Remove the supernatant. Mix the pellet to loosen the cells.



September 2012

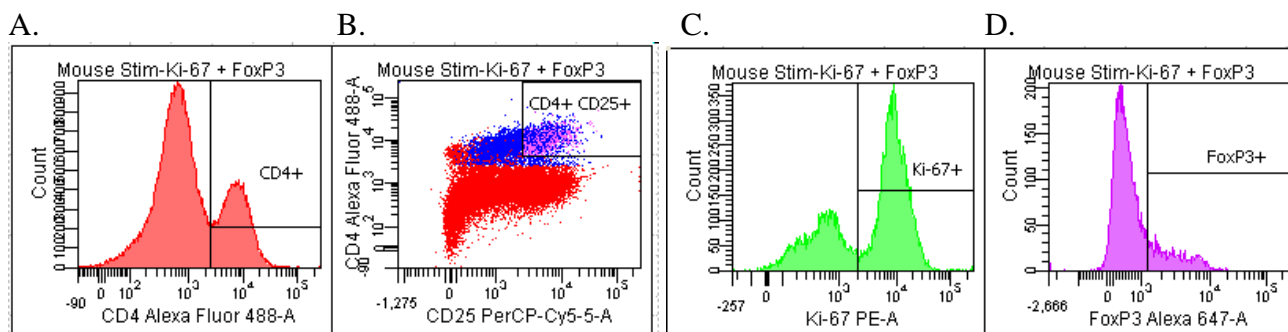
## Simultaneous Measurement of Mouse Foxp3 and Ki-67 in Cultured Splenocytes

### Procedure: Ki-67 Staining

1. Add 250  $\mu$ L of Fix/Perm Solution to each sample tube and mix well. Incubate for 20 minutes at 4°C.
2. Wash the cells twice with 2 mL of 1X Perm/Wash Buffer. Pellet the cells by centrifugation at 500g for 10 minutes.
3. Remove the supernatant, leaving about 100  $\mu$ L in the tube. Mix the pellet to loosen the cells. Note: Perm/Wash Buffer must be maintained in washing steps to keep cells permeabilized.
4. Add 20  $\mu$ L of Anti-Ki-67 antibody to the sample tubes and mix well. Incubate for 30 minutes in the dark at RT.
5. Wash the cells twice with 2 mL of 1X Perm/Wash Buffer. Pellet the cells by centrifugation at 500g for 10 minutes.
6. Remove the supernatant, leaving about 100  $\mu$ L in the tube. Mix the pellet to loosen the cells.
7. Resuspend the cells in 0.5 mL of Stain Buffer and analyze immediately.

### Flow Cytometric Analysis of Stained Cell Samples

Flow cytometric analysis of the samples can be performed by using a flow cytometer equipped with two lasers, a 488-nm blue laser and a 635-nm red laser.

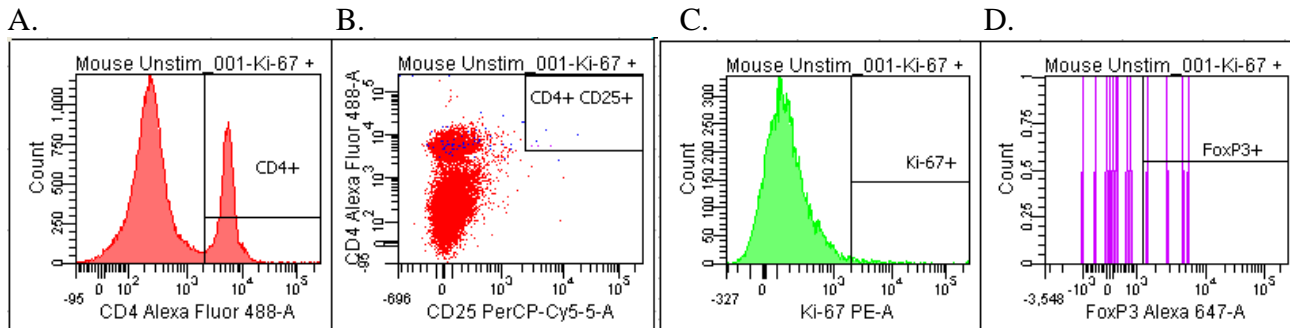


**Figure 1.** Results from the analysis of stimulated cells. The cells were stimulated with anti-CD3 and anti-CD28 antibodies for 72 hours. NA/LE anti-mouse CD3e was coated onto sterile tissue culture plates at a concentration of 10  $\mu$ g/mL in sterile 1X PBS. Soluble NA/LE anti-CD28 was used at 2  $\mu$ g/mL. Panel C shows that 66% of the CD4<sup>+</sup> cells were Ki-67<sup>+</sup>. Panel D shows that 19% of the CD4<sup>+</sup>, CD25<sup>+</sup> cells were Foxp3<sup>+</sup>.

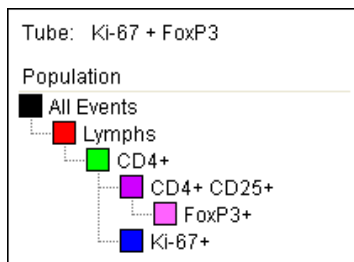


September 2012

## Simultaneous Measurement of Mouse Foxp3 and Ki-67 in Cultured Splenocytes



**Figure 2.** Results from the analysis of unstimulated cells (negative control). Panel C shows an insignificant number of the unstimulated  $CD4^+$  cells were  $Ki-67^+$ . Panel D shows an insignificant number of the unstimulated  $CD4^+$ ,  $CD25^+$  cells were  $Foxp3^+$ .



**Figure 3.** Population hierarchy.

The sample acquisition in this example was done by using the BD LSRFortessa™ flow cytometer, and the analysis using BD FACSDiva™ software.

Alexa Fluor® is a registered trademark of Molecular Probes, Inc.

Cy™ is a trademark of Amersham Biosciences Corp. Cy™ dyes are subject to proprietary rights of Amersham Biosciences Corp and Carnegie Mellon University and made and sold under license from Amersham Biosciences Corp only for research and in vitro diagnostic use. Any other use requires a commercial sublicense from Amersham Biosciences Corp, 800 Centennial Avenue, Piscataway, NJ 08855-1327, USA.

