

## Foxp3 / Transcription Factor Staining Buffer Set

Cat. No. NBP2-266911

Novus' FOXP3 staining kit is optimized for intracellular staining of cells in flow cytometric applications. It is designed and optimized to minimize non-specific staining while maximizing signal-to-noise ratio for clear and consistent data.

### Kit Contents

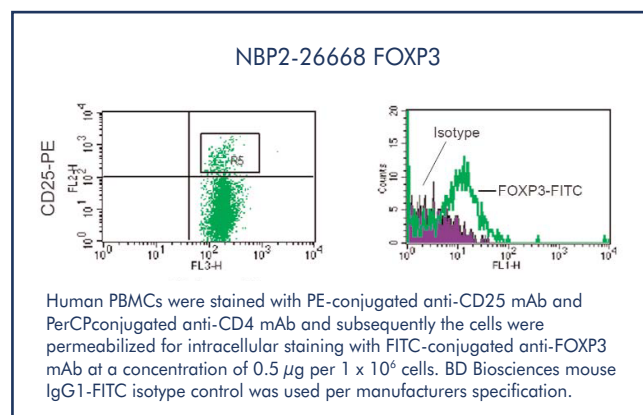
KC-500 Fixation/Permeabilization Concentrate	4X 1 X 15 ml
KC-501 Fixation/Permeabilization Diluent	1X 1 X 50 ml
KC-502 Permeabilization buffer	10X 1 X 50 ml
KC-124 Staining buffer	1X 3 X 60 ml

### Preparation of working solutions

**Fixation/Permeabilization buffer** Prior to staining, dilute the Fixation/Permeabilization Concentrate (1 part) into the Fixation/Permeabilization Diluent (3 parts) to a desired volume of working solution. Working solution should not be stored for more than one day.

**Permeabilization buffer** It is supplied as a 10X solution and can be diluted with deionized water to its final 1X working concentration.

- Determine the number of cells required for staining. For each test sample, it is recommended to use at least  $1 \times 10^6$  cells. The following controls are suggested: a) unstained cells [no primary or secondary antibody staining], b) cells with an isotype control antibody, c) cells with a positive control antibody.
- Harvest the cells and spin down to a pellet at 1000 RPM for 10 min. and decant supernatant.
- Depending on the size of the pellet, resuspend in 2-3 ml of 1X PBS. An exact volume is not necessary at this step.
- Count the cells with a hemocytometer. Remove  $1 \times 10^6$  cells for each sample (including controls) to be tested to a clean conical centrifuge tube. Add 1 ml of 1X PBS to make the decanting easier.
- Centrifuge cells at 1000 RPM for 10 min and decant supernatant.
- Tap the conical tube gently to loosen the pellet.
- Resuspend pellet with the appropriate volume of staining buffer or PBS (50  $\mu$ l per  $1 \times 10^6$  cells).
- Aliquot 50  $\mu$ l of cell suspension to labeled flow tubes.
- Stain surface molecules CD4 and CD25 following the Surface Staining Protocol (Cat. No. NBP2-29481) except for tube with unstained cells. Gently mix and incubate on ice for 30 min in the dark.
- Wash with 2 ml cold Staining buffer (or cold PBS).
- Centrifuge at 1000 RPM for 10 min. and decant supernatant.
- Resuspend cell pellet with pulse vortex and add 1 ml of freshly prepared Fixation/Permeabilization working solution to each sample. Pulse vortex again.
- Incubate at 4°C for 30-60 min in the dark.
- Wash once by adding 2 ml of 1X Permeabilization buffer.
- Centrifuge at 1000 RPM for 10 min. and decant supernatant.
- Optional: Block with 2% mouse serum by adding 2  $\mu$ l to residual buffer after decanting (~100  $\mu$ l) and incubate at 4°C for 15 min.
- Without washing after blocking step, add fluorochrome conjugated anti-FOXP3 antibody or isotype control in 1X Permeabilization buffer (~40  $\mu$ l total) and incubate at 4°C for at least 30 min in the dark. Please perform further titration for optimal staining in your own assay system.
- Wash cells with 2 ml of 1X Permeabilization buffer. Centrifuge and decant supernatant.
- Repeat Step 18.
- Resuspend in appropriate volume Staining buffer and analyze on flow cytometer following manufacturer's recommendations.  
Note: If not analyzing on the same day, samples can be stored overnight, in the dark, at 4°C. Due to the fixation and permeabilization procedure, the FSC/SSC distribution of the cell population will be different than live cells. Therefore the gate and voltages will need to be modified.



Caution: Fixation buffer contains paraformaldehyde which is toxic by inhalation, skin contact, or swallowing. Permeabilization and staining buffers contain 0.05% sodium azide. The 10X Permeabilization buffer has a natural tendency to precipitate, however, its function is not affected by this. To clarify, the solution can be filtered after dilution to 1X working solution. Use caution when handling. All the materials included in this kit should be treated as hazardous materials and be disposed of accordingly.