

Cell Surface Staining Flow Assay Kit

Catalog No. NBP2-26247

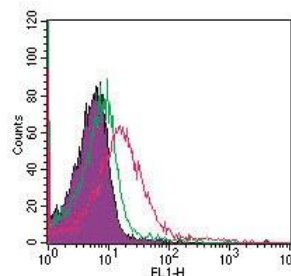
Novus' Cell Surface staining flow kit is a convenient way to quickly and efficiently prepare samples for flow cytometric analysis.

Kit Contents (Store at 4 °C)

KC-136	Staining Buffer	3 x 60 mL
KC-125	Paraformaldehyde 10%	10 mL

Protocol

- Determine the number of cells required for staining. For each test sample, a final concentration of 1×10^6 cells in 50 μ L of staining buffer will be needed. The following controls are suggested:
 - unstained cells [no primary or secondary antibody staining]
 - 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; 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 1×10^6 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.
- Spin down cells to a pellet at 1000 RPM for 10 min and decant supernatant.
- Tap the conical tube gently to loosen the pellet.
- Resuspend pellet with an appropriate volume (50 μ L per 1×10^6 cells) of Staining buffer. Aliquot 50 μ L of cell suspension to individual flow cytometer compatible tubes, one aliquot for each sample to be tested.
- To wash cells, add 1 mL of Staining buffer to each tube, spin down cells to a pellet at 1000 RPM for 10 min and decant supernatant. Prepare primary antibody solutions during centrifugation.
- Dilute each antibody to be tested to the desired concentration in 50 μ L of staining buffer. Resuspend each cell pellet with the appropriate primary antibody. Pipette up and down to thoroughly mix the antibody/cell suspension.
- Incubate on ice for 30 min (protect from light if using a fluorescent labeled primary antibody).
- Centrifuge at 1000 RPM for 10 min and decant supernatant.
- Note: If using a fluorescent-labeled primary antibody, skip Steps 12-14.
- Wash the cells by resuspending each cell pellet with 2 mL of staining buffer, centrifuge at 1000 RPM for 10 min, and decant supernatant. While centrifuging, dilute secondary antibody (FITC, PE or Biotin labeled) in 50 μ L of Staining buffer per sample.
- Resuspend cells with diluted secondary antibody.
- Incubate on ice (protected from light) for 30 min. Centrifuge at 1000 RPM for 10 min and decant supernatant.
- Wash cells twice in 2 mL of Staining Buffer, centrifuging and decanting after each wash step.
- After the final decanting, add 1 mL of Staining Buffer to each tube. Note: If not analyzing on the same day, resuspend cells in 1% paraformaldehyde in Staining Buffer to "fix" cells, and store overnight at 4 °C. The Fixation Buffer can be removed, and the cells prepared for analysis by repeating step 15 and adding 1 mL of Staining buffer to each tube.
- Test samples on a flow cytometer following manufacturer recommendations.



Cell surface flow cytometric analysis of TLR4 in mouse splenocytes using Novus' anti-TLR4 monoclonal antibody (NBP2-24772) and cell surface staining kit. The shaded histogram represents cells alone; green represents isotype control; red represents anti-TLR4 antibody.