

## NBP1-30426 Protocol

### Flow Cytometry protocol for CD3 epsilon Antibody (NBP1-30426)

Flow Cytometry (NBP1-30426): [https://www.novusbio.com/products/cd3-epsilon-antibody-apa1-1\\_nbp1-30426](https://www.novusbio.com/products/cd3-epsilon-antibody-apa1-1_nbp1-30426)

Positive control: Human T cells stimulated with anti-CD3 (MEM-57) antibody (1 mg/ml)

Sample preparation:

At the end of stimulation of T cells, perform staining of surface markers (if required) in PBS + 0.1% BSA for 20 min. on ice.

Wash with PBS and fix with 2% formaldehyde, 30 min on ice.

Wash with PBS and incubate in PBS + 0,1% saponine, 5 min. RT.

Incubate the cells in PBS + 1% BSA + 0.03% saponine, 15 min. on ice.

Incubate with fluorescence-labeled APA1/1 antibody (1-5 mg/ml) in PBS + 1% BSA + 0.03% saponine in dark, 20 min. RT.

Wash with PBS + 1% BSA + 0.03% saponine, resuspend in PBS.