

Data Sheet

Anti-TIGIT Antibody, PE-Labeled

Recombinant, Human Monoclonal

Catalog #: 71228-1

Lot #: 160812 **Conc.:** 0.85 mg/ml

Host Species | Isotype:

Human | IgG1

Formulation: 8 mM Na Phosphate pH 7.4, 110 mM NaCl, 2.2 mM KCl and 20% glycerol.

Stability: Stable for at least 12 months at -80°C. Avoid freeze/thaw cycles. Store diluted antibody with carrier protein (BSA 0.1 – 0.5%) at 4°C, stable for 2 weeks.

References:

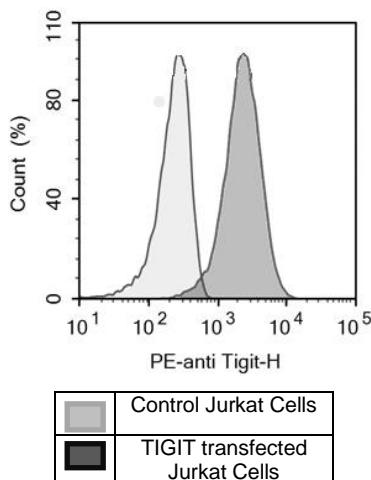
1. Yu, X., *et al.*, *Nat. Immunol.* 2009; **10(1)**: 48-57.
2. Stanitsky, N., *et al.*, *Proc. Natl. Acad. Sci. (USA)*. 2009; **106(42)**: 17858-17863.

Description: R-Phycoerythrin-labeled recombinant human Anti-TIGIT antibody recognizing the TIGIT binding region on immune cells.

Assay Conditions: For flow cytometry, we recommend incubating cells on ice in dark for 15-30 minutes with 20 µg/ml of RPE anti-human TIGIT antibody.

Purification: Protein A affinity chromatography from HEK293 supernatants.

Application: Useful for labeling cells expressing TIGIT for flow cytometry and immunofluorescence microscopy.



Control Jurkat cells and TIGIT-expressing stable Jurkat cells were treated with human PE anti-TIGIT antibody.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**
 Or you can Email us at: info@bpsbioscience.com
 Please visit our website at: www.bpsbioscience.com