

Data Sheet

Anti-TIM-3 Antibody, PE-Labeled

Recombinant, Human Monoclonal
Catalog #: 71221-1
Lot #: 161021 **Conc.:** 2.00

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. Zhu, C. et al., Nat. Immunol. 2005; 6:1245-1252.
2. Anderson, A.C. et al., Science. 2007; 318(5853):1141-1143.

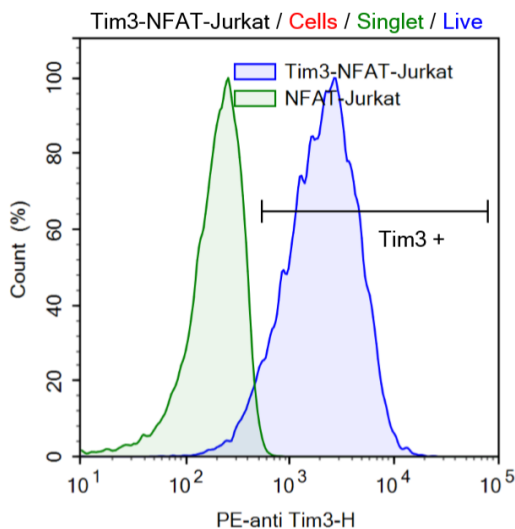
Description: R-Phycoerythrin - labeled recombinant human Anti-TIM-3 antibody recognizing human TIM3 protein. This antibody has not been tested for cross-reactivity with any other species.

Assay Conditions: For flow cytometry, we recommend incubating cells with 20 µg/mL of PE-anti-TIM-3 antibody in 100 µL of FACS buffer for 20 min at 4°C in dark.

Purification: Protein A affinity chromatography from HEK293 supernatants.

Application: Useful for labeling cells expressing TIM-3 for flow cytometry and immunofluorescence microscopy.

Quality Assurance



Control Jurkat cells and TIM-3-expressing stable Jurkat cells were treated with human PE anti-TIM-3 antibody.

Samples	
■	TIM-3-NFAT-Jurkat
■	NFAT-Jurkat

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