

Data Sheet

Anti-Tim3 Antibody

Recombinant, Human Monoclonal

Catalog #: 79098

Lot #: 170718

Conc.: 1.89 mg/ml

Host Species | Isotype:

Human | IgG1

Formulation: 8 mM 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.

References:

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

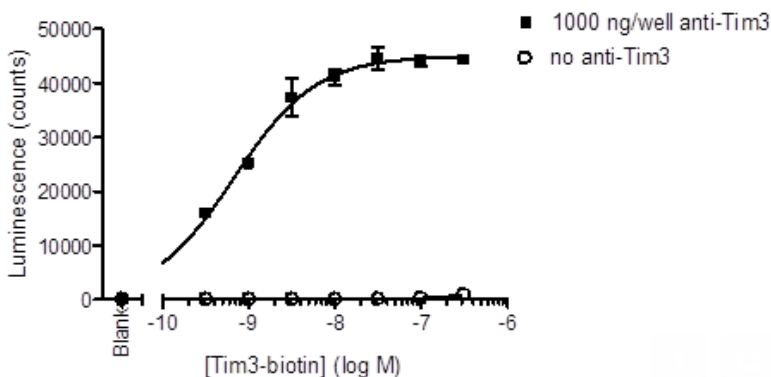
Description: Recombinant human anti-Tim3 antibody recognizing human TIM3 protein. This antibody has not been tested for cross-reactivity with any other species.

Assay Conditions: Tim3-biotin (BPS Bioscience #71237) was titrated from 0 to 300 nM with a 3-fold dilution series in presence of 0 and 1000 ng/well of immobilized anti-Tim3 (BPS Bioscience #79098) and incubated for 1 hour at RT. Binding was detected with Streptavidin-HRP (BPS Bioscience #79742) and ELISA ECL Substrate (BPS Bioscience #79670)

Purification: Protein A affinity chromatography from HEK293 supernatants.

Application: Useful for labeling cells expressing Tim3 for flow cytometry and immunofluorescence microscopy.

Quality Assurance



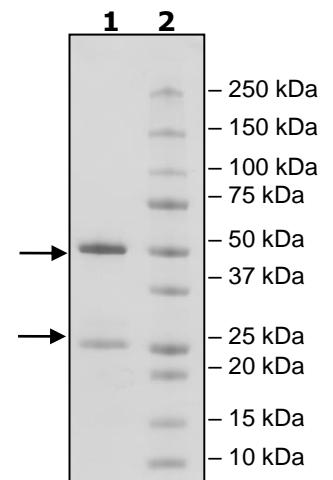
**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
6 µg Anti-Tim3
Antagonist

Lane 2:
Protein Marker

MW:
Heavy chain: 51 kDa
Light chain: 26 kDa

Purity: ≥90%



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

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com