

Data Sheet

TIGIT, Fc fusion, Biotin-labeled

Human, recombinant, Fc fusion protein, with C-terminal AviTag™

Catalog #: 71251

Lot #: 150611

Conc.: 0.79 mg/ml

Formulated in: 8 mM phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol

Stability: At least 6 months at -80°C . Avoid freeze/thaw cycles. Protein may be diluted to $\geq 100 \mu\text{g/ml}$ in PBS + glycerol and stored at -80°C .

References:

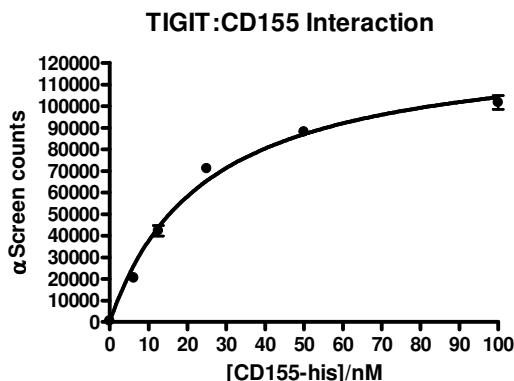
1. Yu, X., *et al.*, *Nat. Immunol.* 2009; **10**(1): 48-57.
2. Stanitsky, N., *et al.*, *Proc. Natl. Acad. Sci. (USA)*. 2009; **106**(42): 17858-17863.

Description: Human T-cell immunoreceptor with Ig and ITIM domains (TIGIT), also known as V-set and immunoglobulin domain-containing protein 9 (VSIG9), and V-set and transmembrane domain-containing protein 3 (VSTM3), GenBank Accession No. NM_173799, a.a. 22-141 fused to Fc region of human IgG1 and enzymatically biotinylated using AviTag™ technology, expressed in a HEK293 cell expression system. MW = 41.6 kDa (monomer). This protein runs at a higher MW by SDS-PAGE due to glycosylation.

Assay Conditions: 100 nM of biotinylated TIGIT was incubated for 1 hour with the indicated concentration of CD155 (Cat. #71181) in a 10 μl reaction in with TIGIT Assay Buffer. Chelate Acceptor beads were added and incubated for 30 min. followed by Strep-donor beads for 30 min and alpha counts were read.

Application: Useful for studying protein binding and for screening small molecules and antibodies.

Quality Assurance

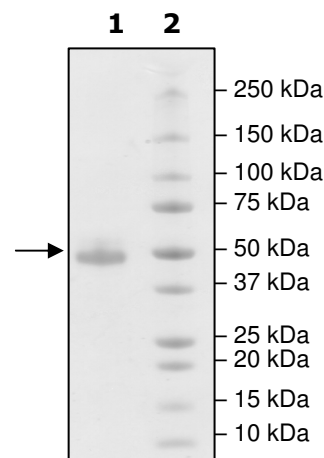


4-20% SDS-PAGE Coomassie staining

Lane 1:
3 μg TIGIT

Lane 2:
Protein Marker

MW: 41.6 kDa
Purity: $\geq 90\%$



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

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

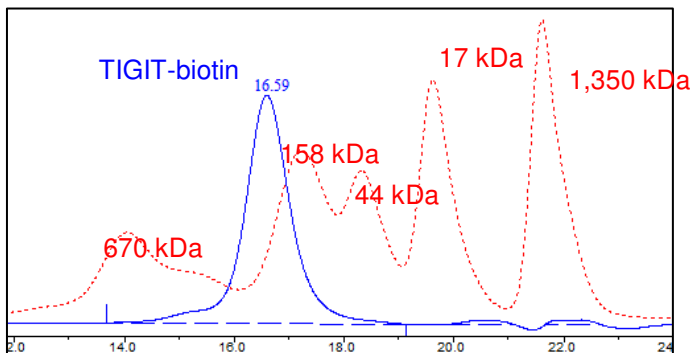
Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6044 Cornerstone Court W, Ste E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Gel Filtration Curve



<10% aggregation

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