

## Data Sheet

### Human CD155, His-tag

Human, recombinant, C-terminal, His-tag

**Catalog #:** 71181

**Lot #:** 160415     **Conc.:** 1.95 mg/ml

**Formulated in:** 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol.

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

**Endotoxin:**  $<1 \text{ EU}/\mu\text{g}$

**References:**

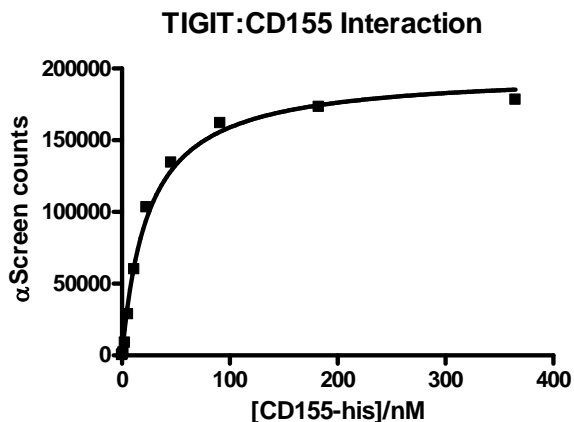
1. Bottino, C. *et al.*, *J. Exp. Med.* 2003; **198(4)**: 557-567.
2. Pende, D. *et al.*, *Blood.* 2005; **105(5)**: 2066-2073.

**Description:** Human CD155, also known as Poliovirus receptor and Nectin-like protein 5, GenBank Accession No. NM\_006505, a.a. 27-343 with a C-terminal His-tag, expressed in a HEK293 cell expression system. MW = 35 kDa. This protein runs at a higher MW by SDS-PAGE due to glycosylation.

**Assay Description:** CD155, His-tag binding to TIGIT was confirmed according to the protocol for BPS Cat. 72029. Briefly, 4 ng of TIGIT-biotin was mixed with varying amounts of CD155, His-tag in TIGIT assay buffer at a final volume of 10  $\mu\text{l}$  and allowed to incubate for 1 hr at room temp. Nickle acceptor beads were added for 30 min, followed by Strep-conjugated acceptor beads for 30 min and reading alpha-counts.

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

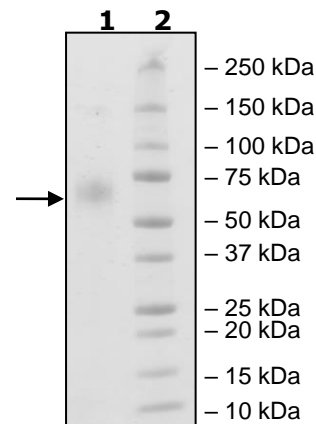
### Quality Assurance



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

**Lane 1:**  
3  $\mu\text{g}$  CD155  
**Lane 2:**  
Protein Marker

**MW:** 35 kDa + glycans  
**Purity:**  $\geq 90\%$



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

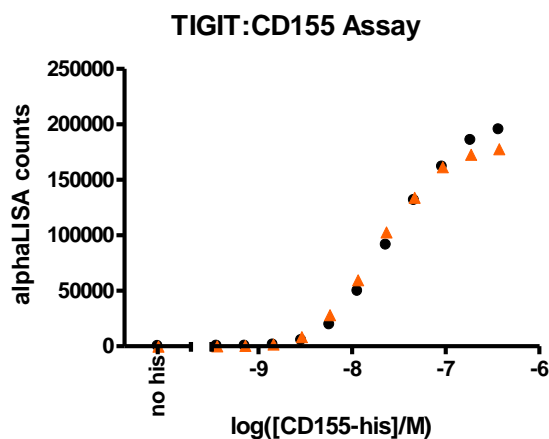
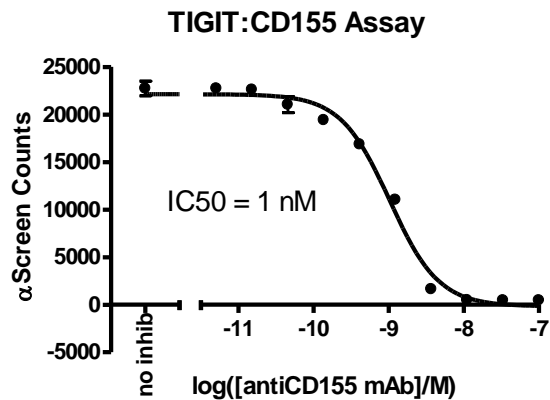
To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**  
 Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)  
 Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)

## Sequence

GDVVVQAPTQVPGFLGDSVTLPCYLQVPNMEVTHVSQLTWARHGEGSGMAV FHQTQ  
 GPSYSESKRLEFVAARLGAELRNASLRMFGLRVEDEGNYTCLFVTFPQGSRSVDIWL  
 RVLAKPQNATAEVQKVQLTGEPVPMARCVSTGGRPPAQITWHS DLGGMPNTSQVPGF  
 LSGTVT V TSLWILVPSSQVDGKNVTCKVEHESFEKPQLLTVNLTVYYPPEVSISGYDN  
 NWYLGQNEATLTCDARSNPEPTGYNWSTTMGPLPPFAVAQGAQLLIRPVDKPINTTLI  
 CNVTNALGARQAELTVQVKEGPPSEHSGMSRNHHHHHHH

Black: CD155 (a.a. 27-343)

Purple: 6x His



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](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)