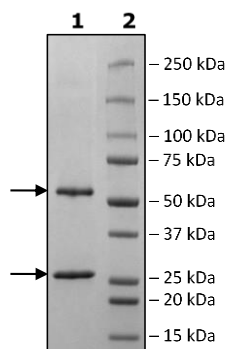


## Product Information

<b>Description:</b>	Human anti-TIGIT neutralizing antibody. This antibody recognizes the region of TIGIT that binds to CD155 and inhibits the interaction between TIGIT and CD155.
<b>Concentration:</b>	3.01 mg/ml
<b>Host Species:</b>	Human
<b>Isotype:</b>	IgG1
<b>Clonality:</b>	Monoclonal
<b>Cross Reactivity:</b>	This antibody has not been tested for cross reactivity with any other species.
<b>Formulated In:</b>	8 mM phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl and 20% glycerol.
<b>Expression System:</b>	HEK293
<b>Format:</b>	Aqueous buffer solution
<b>Stability:</b>	At least 12 months at -80°C. Avoid freeze/thaw cycles.
<b>Storage:</b>	-80°C
<b>MW:</b>	Heavy Chain: 53 kDa, Light Chain: 26 kDa
<b>Purity:</b>	≥90%
<b>Purification:</b>	Protein A affinity chromatography from HEK293 supernatants.
<b>Assay Conditions:</b>	Antibody was added to TIGIT: CD155 homogenous assays at various concentrations. Assay was performed according to the protocol in the TIGIT:CD155 Homogeneous Assay Kit (BPS Bioscience #72029).
<b>Applications:</b>	Useful as a reference antibody for screening TIGIT neutralizing antibodies in biochemical assays.

## Quality Control Data

4-20% SDS-Page Coomassie Staining



TIGIT [Biotinylated]: CD155 Interaction

