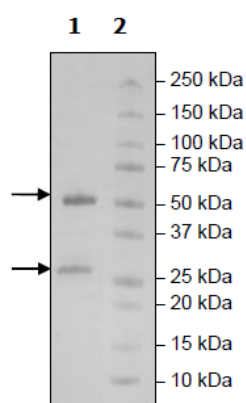


Product Information

Description:	Recombinant human TIGIT inhibitor antibody recognizing an epitope that prevents TIGIT interaction with CD155 on immune cells.
Concentration:	8.23 mg/ml
Host Species:	Human
Clonality:	Monoclonal
Isotype:	IgG1
Formulated In:	8 mM Phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol
Expression System:	HEK293
Format:	Aqueous buffer solution
Stability:	At least 12 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
MW:	Heavy chain: 51 kDa; Light chain: 27 kDa
Purity:	≥90%
Purification:	Protein A affinity chromatography from HEK293 supernatants.
Assay Conditions:	Antibody was added to TIGIT: CD155 homogenous assays at various concentrations. Assay was performed according to the protocol in TIGIT:CD155 Homogeneous Assay Kit (BPS Bioscience #72029)
Applications:	Useful for the inhibition of TIGIT:CD155 and TIGIT: CD112 interactions in biochemical assays. For neutralizing TIGIT in cellular assays, we recommend Anti-TIGIT Neutralizing Antibody (BPS Bioscience #71340).

Quality Control Data

4-20% SDS-Page Coomassie Staining



TIGIT:CD155 Binding Assay

