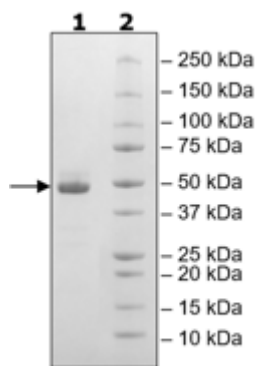


Product Information

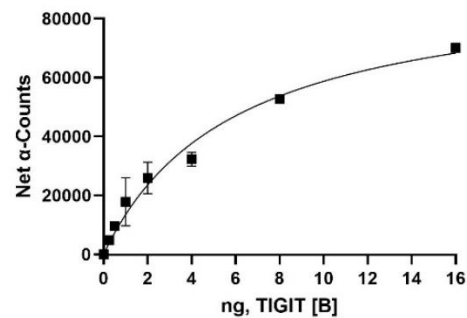
Construct:	TIGIT (22-141-Fc(IgG1)-Avi)-(Biotin)
Label:	This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation is confirmed to be ≥90%.
Concentration:	1.08 mg/ml
Species:	Human
Formulated In:	8 mM phosphate pH 7.4, 110 mM NaCl, 2.2 mM KCl, 20% glycerol
Expression System:	HEK293
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	NM_173799
MW:	42 kDa + glycans
Glycosylation:	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
Purity:	≥90%
Aggregation:	<10%
Assay Conditions:	Assay was performed according to the protocol of the TIGIT:CD155 Homogeneous Assay Kit (BPS Bioscience #72029) with TIGIT titrated from 16 ng/rxn to 0 ng/rxn.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

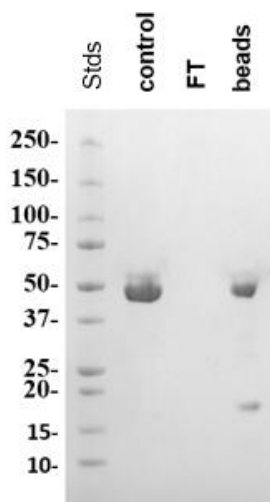
4-20% SDS-Page Coomassie Staining



TIGIT [Biotinylated]: CD155 Interaction



Biotin-Avidin Pulldown



Gel Filtration Trace

