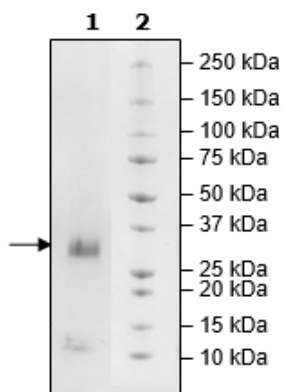


Product Information

Construct:	CD70 (His-Avi-39-193(end))
Concentration:	0.47 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_001252
MW:	22 kDa + glycans
Glycosylation:	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
Purity:	85%
Assay Conditions:	Assay was performed according to the CD27: CD70[Biotinylated] Inhibitor Screening Assay Kit (BPS Bioscience #79695) with CD27 [Biotinylated] enzyme titrated from 300 ng/rxn – 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



CD70: CD27[Biotinylated] Interaction

