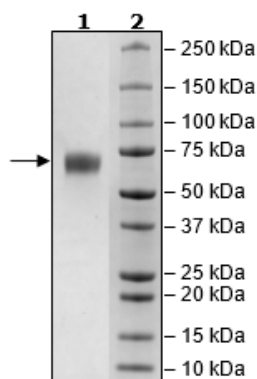


Product Information

Construct:	CD28 (19-152-Fc(IgG1)-Avi)
Concentration:	2.29 mg/ml
Species:	Human
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 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	NM_006139
MW:	41 kDa + glycans
Glycosylation:	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
Purity:	≥90%
Assay Conditions:	Assay was performed according to CD28:B7-1[Biotinylated] Inhibitor Screening Assay Kit (BPS Bioscience #72007) with various amounts of B7-1 [Biotinylated].
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



CD28:B7-1[Biotinylated] Activity

