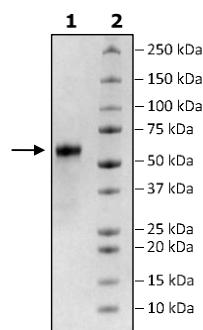


Product Information

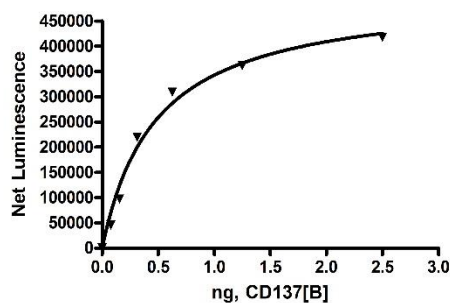
Construct:	CD137 (4-1BB) (24-186-Fc(IgG1)-Avi)-(biotin)
Label:	This protein is enzymatically biotin-labeled using Avi-tag™ technology. Biotinylation is confirmed to be ≥90%.
Concentration:	1.52 mg/ml
Species:	Human
Formulated In:	8 mM phosphate pH 7.4, 110 mM NaCl, 2.2 mM KCl, 20% glycerol
Expression System:	HEK293
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	NM_001561
MW:	46 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 CD137[Biotinylated]:CD137L Inhibitor Screening Assay Kit (BPS Bioscience, #72025) with CD137 [Biotinylated] enzyme titrated from 80 ng/rxn – 0.08 ng/rxn.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

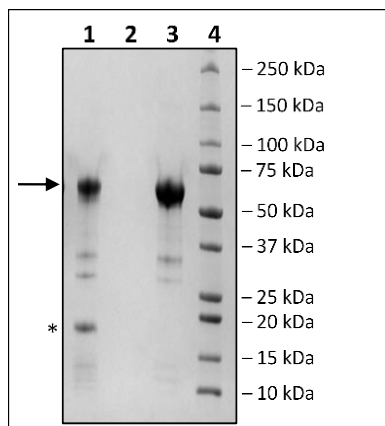
4-20% SDS-Page Coomassie Staining



CD137[Biotinylated]:CD137L Interaction



Biotin-Avidin Pulldown



Lane 1: Beads
Lane 2: FT
Lane 3: Control
Lane 4: Standards

Gel Filtration

