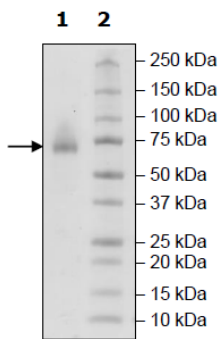


Product Information

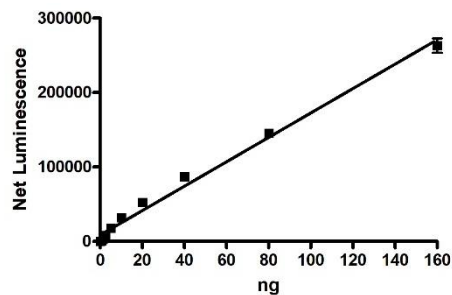
Construct:	CD48, (27-220-Fc(IgG1)-Avi)-(Biotin)
Concentration:	3.02 mg/ml
Species:	Human
Formulated In:	8 mM Phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol.
Expression System:	HEK293
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	NM_001778
MW:	51 kDa +glycans
Purity:	≥90%
Assay Conditions:	Assay was performed according to the CD137[Biotinylated]:CD137L Inhibitor Screening Assay Kit (BPS Bioscience, #72025). The plate was coated with 100 ng of CD244. CD48, Biotin-labeled was titrated from 160 ng/rxn – 0 ng/rxn.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.
Notes:	This protein runs at a higher MW by SDS-PAGE due to glycosylation.

Quality Control Data

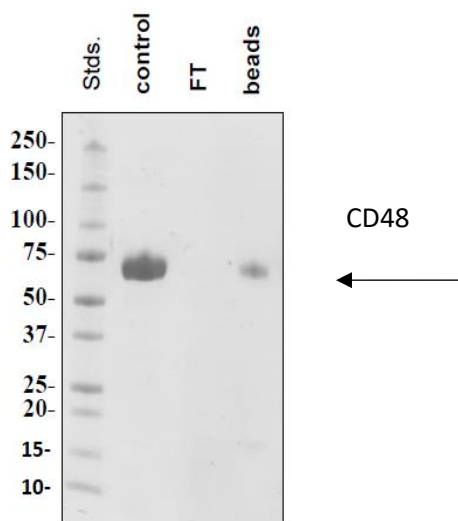
4-20% SDS-Page Coomassie Staining



CD244: CD48, Biotin-Labeled Activity



Biotin-Avidin Pulldown



This protein is enzymatically biotinylated using Avi-tag™ technology. Biotinylation is confirmed to be $\geq 90\%$.

Gel Filtration Trace

<10% Aggregation

