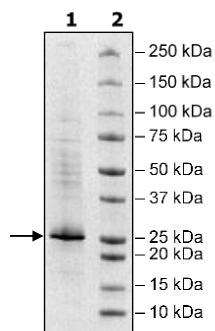


Product Information

Construct:	CD137L (His-50-254(end))
Concentration:	0.91 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_003811
MW:	24 kDa
Purity:	74%
Assay Conditions:	Assay was done according to the CD137[Biotinylated]:CD137L Inhibitor Screening Assay Kit (BPS Bioscience, #72025) with CD137[Biotinylated] titrated from 80 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



CD137[Biotinylated]:CD137L Interaction

