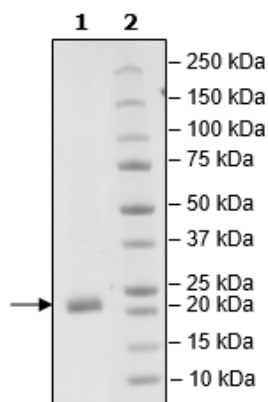


Product Information

Construct:	LIGHT (His-74-240(end))
Concentration:	0.68 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_003807
MW:	21 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 LIGHT:HVEM [Biotinylated] Inhibitor Screening Assay Kit (BPS Bioscience #79684) protocol. Coat protein (LIGHT) was added ON at 4 deg 100 ng/well. Various concentrations of biotinylated ligand (HVEM[B], 0-80 ng) added for 2h at RT. Strep-HRP added for 1h at RT.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-Page Coomassie Staining



LIGHT: HVEM [Biotinylated] Binding

