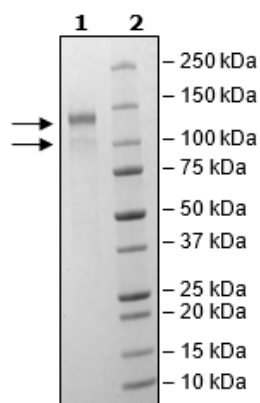


## Product Information

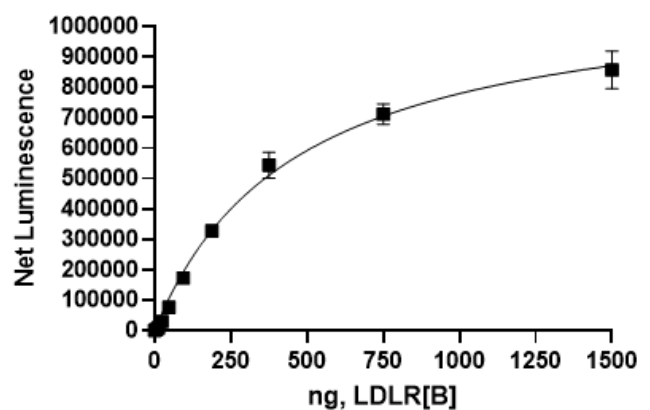
<b>Description:</b>	Recombinant human LDLR (Low Density Lipoprotein Receptor), encompassing amino acids 22-788. This construct contains a C-terminal FLAG-tag followed by an Avi-Tag™. The recombinant protein was enzymatically biotinylated using the Avi-Tag™ and affinity purified.
<b>Species:</b>	Human
<b>Construct:</b>	LDLR (22-788(end)-FLAG-Avi)-(Biotin)
<b>Concentration:</b>	0.36 mg/ml
<b>Expression System:</b>	HEK293
<b>Purity:</b>	≥90%
<b>Format:</b>	Aqueous buffer solution.
<b>Formulated In:</b>	8 mM phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol
<b>MW:</b>	88 kDa + glycans
<b>Glycosylation:</b>	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
<b>Genbank Accession:</b>	NM_000527
<b>Label:</b>	This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation is confirmed to be ≥90%
<b>Stability:</b>	At least 6 months at -80°C.
<b>Storage:</b>	-80°C
<b>Instructions for Use:</b>	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.
<b>Assay Conditions:</b>	Assay was performed according to PCSK9 [Biotinylated]-LDLR Binding Assay Kit (BPS Bioscience #72002) with PCSK9 coated at 4 ng/μl and LDLR[Biotinylated] titrated at various concentrations.
<b>Applications:</b>	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

## Quality Control Data

4-20% SDS-PAGE Coomassie Staining



PCSK9-LDLR[Biotinylated] Interaction



## Quality Control Data

### Biotin-Avidin Pulldown

