

PCSK9, C-terminal His-Avi-Tag, Biotin-Labeled Recombinant

Catalog: 71304

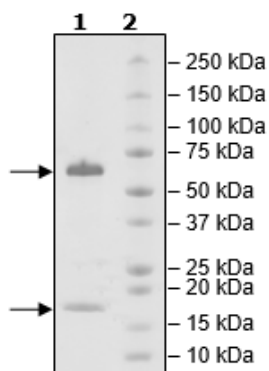
Lot: 211118

Product Information

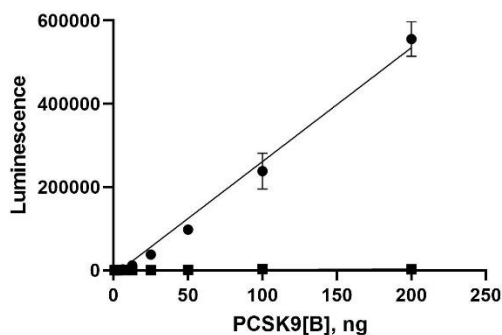
Construct:	PCSK9 (31-692(end)-His-Avi)-(Biotin)
Label:	This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation confirmed to be ≥90%.
Concentration:	1.67 mg/ml
Species:	Human
Formulated In:	8 mM phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol
Expression System:	HEK293
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	Pro-PCSK9: PCSK9: NM_174936
MW:	74 kDa
Purity:	≥90%
Agregation:	<10%
Assay Conditions:	Assay was performed according to PCSK9[Biotinylated]-LDLR Binding Assay Kit (BPS Bioscience #72002), with the PCSK9[Biotinylated] titrated from 200 ng- 0 ng.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

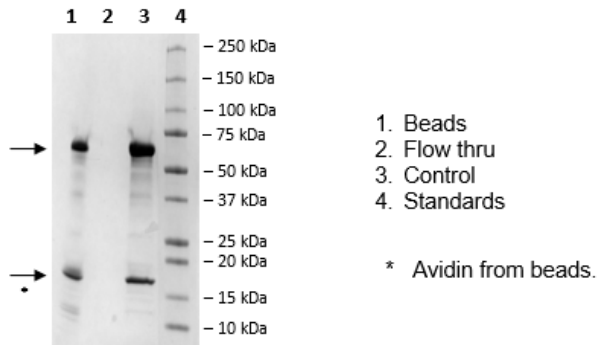
4-20% SDS-Page Coomassie Staining



PCSK9[Biotinylated]:LDLR Activity



Biotin-Avidin Pulldown



Gel Filtration Trace

