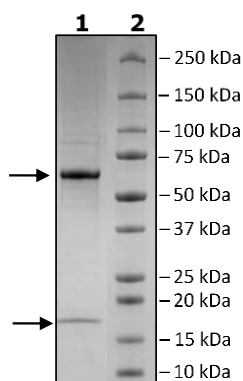


Product Information

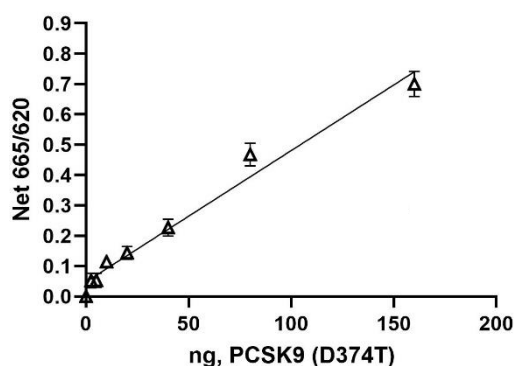
Construct:	PCSK9 (D374T) (31-692(end)-His-Avi)-(Biotin)
Mutation:	D374T
Label:	This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation confirmed to be ≥90%.
Concentration:	0.90 mg/ml
Species:	Human
Formulated In:	8 mM phosphate pH 7.4, 110 mM NaCl, 2.2 mM KCl, 20% glycerol
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_174936
MW:	Pro: 14 kDa + glycans; Mature: 60 kDa + glycans
Glycosylation:	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
Purity:	≥90%
Aggregation:	<10%
Assay Conditions:	Assay was done according to the PCSK9(D374T)-LDLR TR-FRET Assay Kit (BPS Bioscience #72011) with PCSK9 (D374T) titrated from 160 ng-2.5 ng/rxn.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

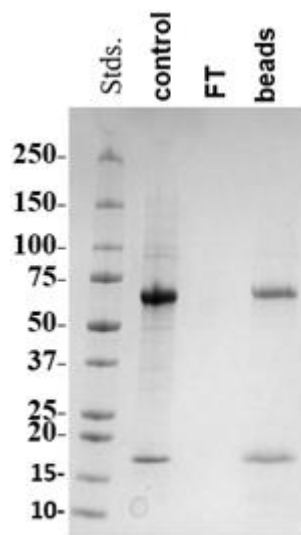
4-20% SDS-Page Coomassie Staining



PCSK9 (D374T): LDLR Activity



Biotin-Avidin Pulldown



Gel Filtration Trace

