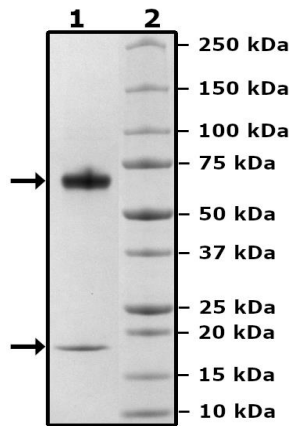


Product Information

Description:	Recombinant human PCSK9 (Human proprotein convertase subtilisin/kexin 9) encompassing amino acids 31-692, which comprise the protodomain, the catalytic, and the C-terminal domains. The construct contains a C-terminal His-tag (6xHis) followed by an Avi-tag™. The protein contains the potent gain of function D374Y mutation. The protein was affinity purified.
Background:	PCSK9 (Proprotein convertase subtilisin/kexin type 9) functions as a negative regulator of hepatic low-density lipoprotein receptors (LDLRs) by binding to the LDLR ectodomain. The D374Y mutation is a gain of function mutation. The D374Y mutation is associated with hypercholesterolemia. This PCSK9 mutant is more potent at decreasing LDL uptake than wild-type PCSK9, most likely by increasing the binding affinity of PCSK9 for the LDLR. A deeper understanding of the function of this mutant protein and development of specific inhibitors may result in new therapeutic approaches for PCSK9 D374Y-linked diseases.
Species:	Human
Construct:	PCSK9 (D374Y) (31-692(end)-His-Avi)(Biotin)
Mutation:	D374Y
Concentration:	0.44 mg/ml
Expression System:	HEK293
Purity:	≥90%
Format:	Aqueous buffer solution.
Formulated In:	8 mM phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol
MW:	Pro: 14 kDa; Mature: 60 kDa + glycans
Glycosylation:	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
Genbank Accession:	NM_174936
Label:	This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation was confirmed to be ≥90%.
Stability:	At least 6 months at -80°C.
Storage:	-80°C
Instructions for Use:	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.
Applications:	SDS-PAGE and pull-down assays.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



Biotin-Avidin Pulldown

