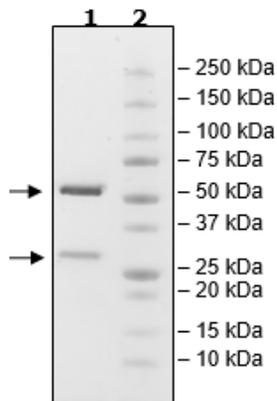


Product Information

Description:	This anti-PCSK9 (proprotein convertase subtilisin/kexin type 9) antibody is a purified, recombinant monoclonal antibody that recognizes human PCSK9. This antibody has been tested for neutralization of PCSK9 interaction with LDLR (low density lipoprotein receptor) in a binding assay. The heavy chain of this antibody contains a C-terminal Avi-Tag™.
Background:	PCSK9 (Proprotein convertase subtilisin/kexin type 9) functions as a negative regulator of hepatic low-density lipoprotein receptors (LDLRs) and therefore is a critical regulator of cholesterol metabolism. It is an endopeptidase that binds to the EGFR-like ectodomain of LDLR, leading to LDLR degradation, which in turn, results in increased circulating LDL. Inhibiting the PCSK9-LDLR interaction is an increasingly desirable therapeutic approach for lowering LDL-cholesterol levels to replace or supplement statins. New therapies are critical for addressing atherosclerosis, stroke, heart disease, and other cardiovascular disorders.
Species:	Human
Isotype:	IgG1
Clonality:	Monoclonal
Concentration:	0.22 mg/ml
Expression System:	HEK293
Purification:	Protein A affinity chromatography from HEK293 supernatants.
Purity:	≥90%
Format:	Aqueous buffer solution.
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol
MW:	Heavy Chain: 54 kDa; Light Chain: 25 kDa
Stability:	At least 12 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.
Assay Conditions:	Assay was performed according to the PCSK9-LDLR Homogeneous Assay Kit (BPS Bioscience #78812) with the Anti-PCSK9 Neutralizing Antibody, Biotin-labeled diluted three-fold from 200 nM to 0 nM.
Applications:	Useful for the studying the binding and neutralization of PCSK9 to LDR in ELISA and cellular assays.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



PCSK9:LDLR Binding

