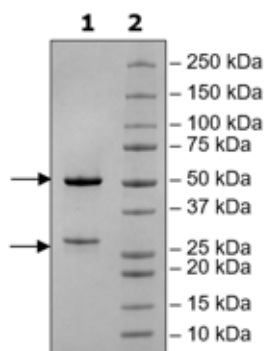


Product Information

Description:	Neutralizing recombinant human antibody (IgG1) recognizing the LDLR binding region of human PCSK9.
Concentration:	2.32 mg/ml
Species:	Human
Clonality:	Monoclonal
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
MW:	Heavy Chain: 50 kDa; Light Chain: 23 kDa
Purity:	≥90%
Purification:	Protein A affinity
Assay Conditions:	Assay was performed according to the PCSK9[Biotinylated]-LDLR Binding Assay Kit (BPS Bioscience #72002) with PCSK9 [Biotinylated] enzyme concentrated at 2.5 ng/μl and the Anti-PCSK9 neutralizing antibody diluted three-fold from 100 nM to 0 nM.
Applications:	Useful as an inhibitor of PCSK9-LDLR binding using the PCSK9[Biotinylated]-LDLR Binding Assay Kit (BPS Bioscience #72002).

Quality Control Data

4-20% SDS-Page Coomassie Staining



PCSK9 [Biotinylated]: LDLR Binding Interaction

