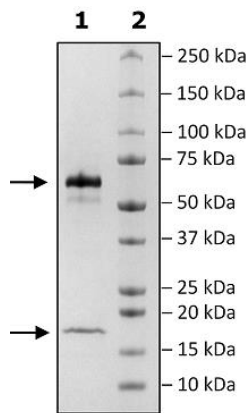


Product Information

| | |
|---------------------------|--|
| Construct: | PCSK9(31-692(end)-His) |
| Concentration: | 1.02 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; Mature: 58 kDa + glycans |
| Glycosylation: | This protein runs at higher MW by SDS-PAGE due to glycosylation. |
| Purity: | ≥90% |
| Assay Conditions: | Assay was performed according to the PCSK9[Biotinylated]-LDLR Binding Assay Kit (BPS Bioscience, #72002) with PCSK9 coated instead and LDLR[Biotinylated] enzyme titrated instead of LDLR from 400 ng/rxn – 0.39 ng/rxn. |
| Applications: | Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling. |
| Notes: | PCSK9 is autocleaved to the ~14 kDa prodomain (a.a. 31-152) and the ~60 kDa mature form (a.a. 153-692(end)), which run at higher apparent MW by SDS-PAGE. |

Quality Control Data

4-20% SDS-Page Coomassie Staining



PCSK9:LDLR[Biotinylated] Binding

