

Data Sheet

PCSK9, Biotinylated, His-tags

Human, recombinant, C-terminal His-Avi Tags

Catalog #: 71304

Lot #: 150318

Conc.: 0.99 mg/ml

Formulated in: 8 mM Phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol

Stability: At least 6 months at -80°C . Avoid multiple freeze/thaw cycles. Storing diluted protein is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).

References:

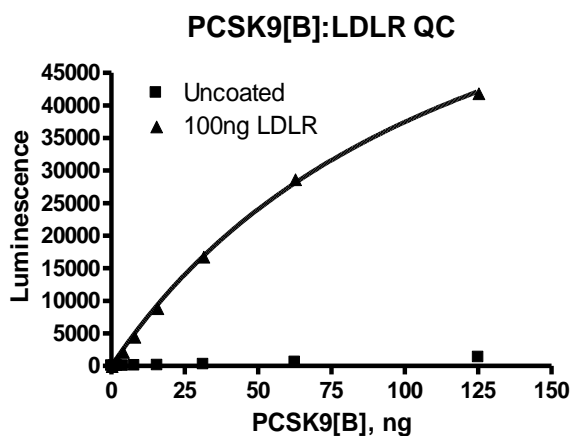
1. Abifadel, M., *et al.*, *Nat Genet.* 2003; **34(2)**: 154-6.
2. Abifadel, M., *et al.*, *Hum Mutat.* 2009; **30(4)**: 520-9.

Description: Human proprotein convertase subtilisin/kexin type 9 (PCSK9), also known as FH3, HCHOLA3, and PC9, GenBank Accession No. NM_174936, a.a. 31-692(end), with C-terminal His and Avi tags, MW=73.8 kDa (calculated), expressed in an HEK293 cell expression system and enzymatically biotinylated using Avitag™ technology. PCSK9 is autocleaved to the ~14 kDa pro-domain (a.a. 31-152) and the ~60 kDa mature form (a.a. 153-692(end)), which run at higher apparent MW by SDS-PAGE.

Assay Conditions: 50 mL reaction mix (50 mM HEPES pH 7.4, 50 mM NaCl, 300 nM CaCl_2 , 0.05% Tween, 0.1% BSA, and 0–1500 ng Biotin-Labeled PCSK9) is added to wells coated with LDLR ectodomain. Incubate for 2 hr. Add streptavidin-labeled HRP and incubate 1 hr. Finally, add HRP chemiluminescent substrates and read luminescence.

Application: Useful for studying protein binding and screening small molecules.

Quality Assurance

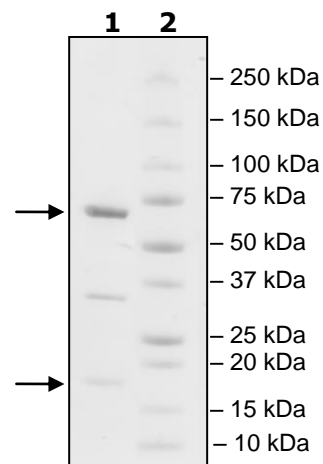


**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
4 µg PCSK9

Lane 2:
Protein Marker

MW: 73.8 kDa
Purity: ≥80%



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6044 Cornerstone Court W, Ste E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Sequence:

QEDEDGDYEELVLALRSEEDGLAEAPEHGTTATFHRCADPWRLPGTYVVVLKEETHL
SQSERTARRLQAQAARRGYLTKILHVFHGLLPGLVKMSGDLLELALKLPHVDYIEEDS
SVFAQSIPWNLERITPPRYRADEYQPPDGGSLVEVYLLDTSIQSDHREIEGRVMVTD
NVPEEDGTRFHRQASKCDSHGTHLAGVVSGRDAGVAKGASMRSLRVLNCQGKGT
SGTLIGLEFIRKSQLVQPVGPLVLLPLAGGYSRVLNAACQRLARAGVVLVTAAGNFR
DDACLYSPASAPEVITVGATNAQDQPVTLGTLGTNFGRCVDLFAPGEDIIIGASSDCST
CFVSQSGTSQAAAHVAGIAAMMLSAEPELTLAELRQRLIHFSKDVINEAWFPEDQR
VLTPNLVAALPPSTHGAGWQLFCRTVWSAHS GPTRMATAVARCAPDEELLSCSSFS
RSGKRRGERMEAQGGKLCRAHNAFGGEGVYAIARCCLLPQANCSVHTAPPAEAS
MGTRVHCHQQGHVLTGCSSHWEVEDLGTHKPPVLRPRGQPNQCVGHREASIHASC
CHAPGLECKVKEHGIPAPQEQT VACEEGWTLTGCSALPGTSHVLGAYAVDNTCVV
RSRDVSTTGSTSEGAVTAVAICCRSRHLAQASQELQH H H H H H H G G G L N D I F E A Q K I E W
HE

Green: Pro-PCSK9, a.a. 31-152
Black: PCSK9, a.a. 153-692(end)
Red: His tag
Blue: Avi Tag

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com