

Data Sheet

PCSK9, Biotinylated

Human, recombinant, C-terminal Avi-His Tags

Catalog #: 71305

Lot #: 141021-BD1

Conc.: 1.77 mg/ml

Formulated in: 40 mM Tris, pH 8.0, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol.

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

References:

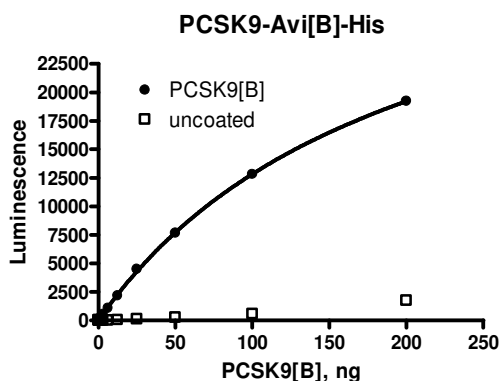
1. Abifadel, M., *et al.*, *Nat Genet.* 2003 Jun;**34**(2):154-6.
2. Abifadel, M., *et al.*, *Hum Mutat.* 2009 Apr;**30**(4):520-9.

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 Avi and His 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, 200 μM CaCl_2 , 0.05% Tween, 0.1% BSA, and 0–200 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

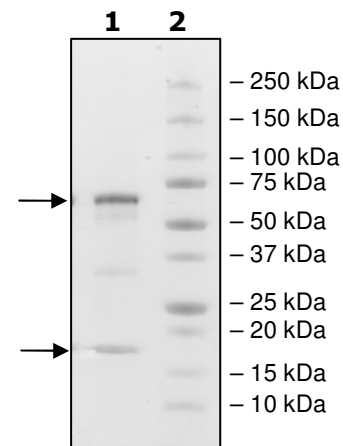


**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
4 μg PCSK9

Lane 2:
Protein Marker

MW: 73.8 kDa
Purity: $\geq 80\%$



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

Protein Sequence:

QEDEDGDYEELVLALRSEEDGLAEAPEHGTTATFHRCADPWRLPGTYVVVLKEETHL
SQSERTARRLQAQAARRGYLTKILHVFHGLLPGLVKMSGDLLELALKLPHVDYIEEDS
SVFAQSIPWNLERITPPRYRADEYQPPDGGSLVEVYLLDTSIQSDHREIEGRVMVTD
NVPEEDGTRFHRQASKCD SHGTHLAGVVSGRDAGVAKGASMRSLRVLNQCQKGTVS
GTLIGLEFIRKSQLVQPVGPLVVLLPLAGGYSRVLNAACQRLARAGVVLVTAAGNFRDD
ACLYSPASAPEVITVGATNAQDQPVTLGTLGTNFGRCVDFAPGEDIIIGASSDCSTCFV
SQSGTSQAAAHVAGIAAMMLSAEPELTLAELRQRLIHFSKDVINEAWFPEDQRVLT
LVAALPPSTHGAGWQLFCRTVWSAHS GPTRMATAVARCAPDEELLSCSSFSRSGKRR
GERMEAQGGKLCRAHNAFGGEGVYAIARCCLLPQANCSVHTAPPAEASMGTRVHC
HQQGHVLTGCSSHWEVEDLGTHKPPVLRPRGQPNQCVGHREASIHASCCHAPGLEC
KVKEHGIPAPQEQVTVACEEGWTLTGCSALPGTSHVLGAYAVDNTCVVRSRDVSTTG
STSEGAVTAVAICCRSRHLAQASQELQGGGLNDIFEAQKIEWHEHHHHH

Green: Pro-PCSK9 31-152

Black: PCSK9 a.a. 153-692(end)

Blue: AviTag

Red: His 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