

Data Sheet

PCSK9, His-Avi-Tag, Biotin-Labeled

Human, Recombinant, C-terminal His-Avi Tags

Catalog #: 71304-2

Lot #: 170126-3 **Conc.:** 0.76 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.

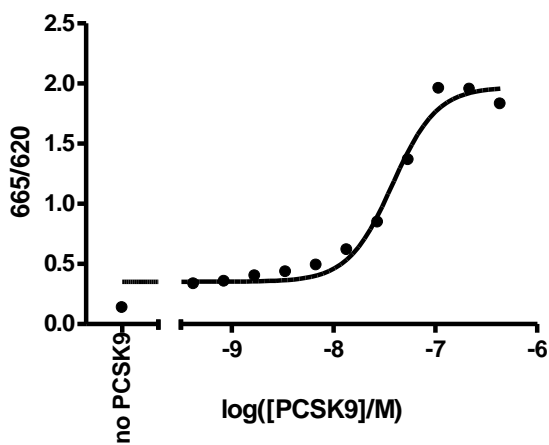
Description: Human proprotein convertase subtilisin/kexin type 9 (PCSK9), also known as FH3, HCHOLA3, and PC9, GenBank Accession No. NM_174936, a.a. 31-692(end), with C-terminal His and Avi tags, MW=74 kDa (calculated), expressed in an HEK293 cell expression system and enzymatically biotinylated using Avitag™ technology. 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.

Assay Conditions: Mixed Streptavidin-XL (1x) with LDLR-Eu (1.5 nM) and initiated the binding reaction by adding various concentrations of PCSK9 biotin (0-430 nM respectively). Incubated for 2h and measured TR-FRET 665/620.

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

Quality Assurance

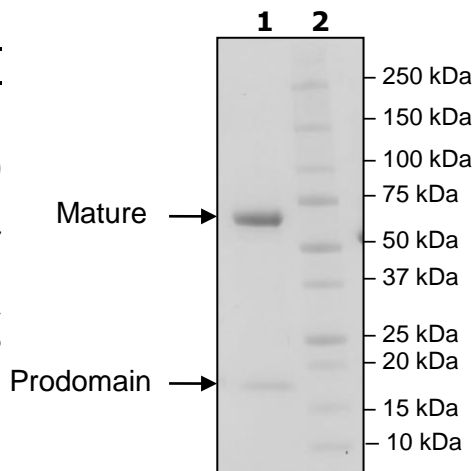
Specific Activity



4-20% SDS-PAGE
Coomassie staining

Lane 1: 5µg PCSK9
Lane 2: Protein Marker

MW: 74 kDa
Purity: ≥90%



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**

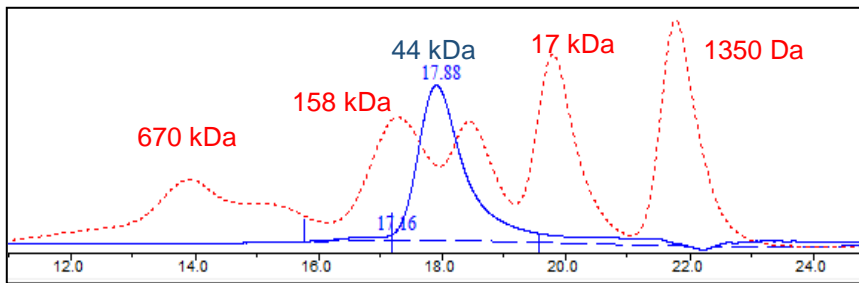
Or you can Email us at: info@bpsbioscience.com

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

Gel Filtration Trace



<10% Aggregation

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