

Data Sheet

PCSK9, His-Avi-Tag (Mouse)

Mouse, Recombinant, C-terminal
His-Avi-tag

Catalog #: 71221-1

Lot #: 130110-D **Conc.:** 0.24 mg/ml

Formulated in: 40 mM Tris-HCl pH 8.0,
110 mM NaCl, 2.2 mM KCl, 0.04%
Tween20, 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: Mouse proprotein convertase subtilisin/kexin type 9 (PCSK9), also known as FH3, HCHOLA3, and PC9, GenBank Accession No. NM_153565, a.a. 35-694 (end), with C-terminal His-Avi-tag, MW=74 kDa (calculated), expressed in a HEK293 cell expression system.

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

Quality Assurance

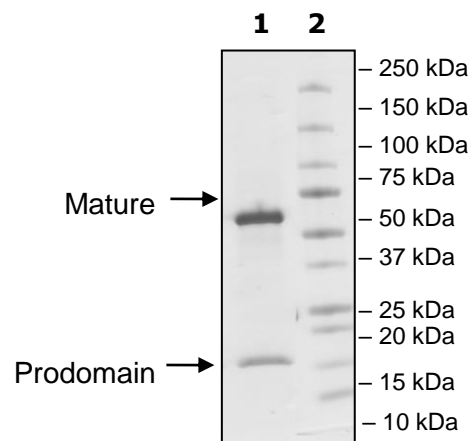
4-20% SDS-PAGE Coomassie staining

Lane 1:
4 μg PCSK9

Lane 2:
Protein Marker

MW: 74 kDa

Purity: $\geq 77\%$



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