

Data Sheet

PCSK9, His-Avi-Tag (Rat)

Rat, Recombinant, C-terminal His- tag
Catalog #: 71220-2
Lot #: 130118-Dc1 **Conc.:** 0.39 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: Rat proprotein convertase subtilisin/kexin type 9 (PCSK9), also known as FH3, HCHOLA3, and PC9, GenBank Accession No. NM_199253, a.a. 31-691(end), with C-terminal His-Avi-tag, MW=76 kDa (calculated), expressed in a HEK293 cell expression system.

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

Quality Assurance

1

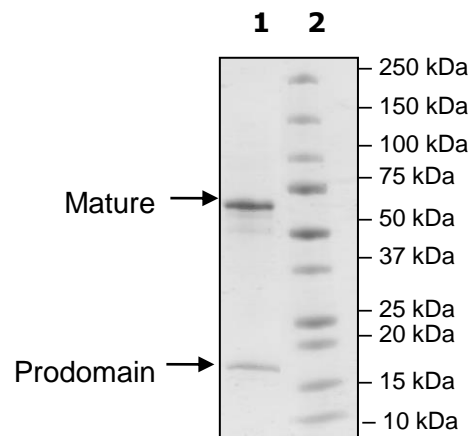
**4-20% SDS-PAGE
 Coomassie staining**

Lane 1:
 4 μg PCSK9

Lane 2:
 Protein Marker

MW: 76 kDa

Purity: $\geq 67\%$



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