

## Data Sheet

### PCSK9(D374T), Biotinylated

Human, recombinant, C-Terminal His-Avi Tags

**Catalog #:** 71211

**Lot #:** 150203-D

**Conc.:** 0.68 mg/ml

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

**Stability:** At least 6 months at  $-80^{\circ}\text{C}$ . Avoid freeze/thaw cycles. Storing diluted enzyme is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).

**References:**

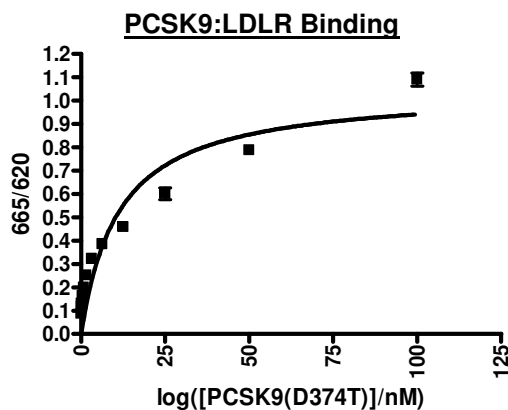
1. Chan, J.C. *et al.* (2009). *Proc. Natl. Acad. Sci. USA*, **106**: 9820-9825.
2. Liang, H., *et al.* (2012). *J. Pharmacol. Exp. Ther.* **340**: 2289-236.

**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 and a D374T mutation, MW=73.8 kDa, 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), which run at higher MW by SDS-PAGE.

**Assay Conditions:** Reaction was performed using PCSK9(D374T)-LDLR TR-FRET Assay Kit (Cat. #72011). Briefly, 5 ng of europium-labeled LDLR were mixed with varying amounts of PCSK9(D374T) and dye-labeled acceptor in PCSK9 TR-FRET assay buffer and incubated at RT for 2 hr. Donor emission was read at 620 nm followed by acceptor emission at 665 nm.

**Application:** Useful for the study of enzyme kinetics, binding studies, screening inhibitors, and selectivity profiling.

### Quality Assurance



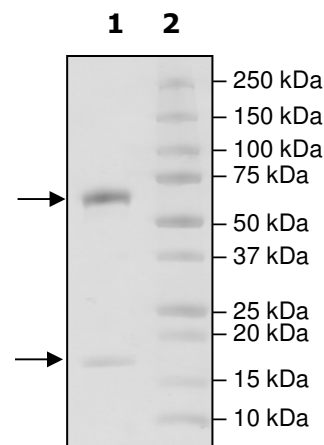
**4-20% SDS-PAGE  
Coomassie staining**

**Lane 1:**  
3 µg PCSK9

**Lane 2:**  
Protein Marker

**MW:** Pro: 14 kDa  
Mature: 60 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](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)