



6044 Cornerstone Court W, Ste E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Data Sheet

Anti-PCSK9, Biotin-labeled

Recombinant, Human Monoclonal
Anti-PCSK9

Catalog #: 71216

Lot #: 141027-BDS **Conc.:** 0.22 mg/ml

Host Species | Isotype:

Human | IgG1

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

Endotoxin Level: <1 EU/μg

Stability: Stable for at least 12 months at
-80°C. *Avoid freeze/thaw cycles. Store
diluted antibody with carrier protein (BSA 0.1
– 0.5%) at 4°C, stable for 2 weeks.*

References:

1. Banerjee, Y., *et al.*, *N. Engl. J. Med.*
2012; **366(25)**:2425-2426.
2. Maxwell, K.N., *Circ. Res.* 2012; **111(3)**:
274-277.

Description: Purified recombinant human
anti-PCSK9 (IgG1) antibody with a C-terminal
Avi-tag, expressed in HEK293 cells. Purified
antibody was enzymatically biotinylated using
Avitag™ technology. This antibody reacts with
human PCSK9 and has not been tested for
cross-reactivity with other species.

Purification: Protein A affinity
chromatography from HEK293 supernatants.

Applications: Western blot, ELISA

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