

# Product Information



## PCSK9 (mouse) Polyclonal Antibody

Item No. 10008811 • Lot No. XXXXX

<b>Contents:</b>	This vial contains <i>lot specific</i> mg of peptide affinity-purified IgG in <i>lot specific</i> µl TBS, pH 7.4, 50% glycerol, 0.5 mg/ml BSA, 0.02% sodium azide.
<b>Synonyms:</b>	NARC-1; Proprotein Convertase Subtilisin Kexin 9
<b>Antigen:</b>	Synthetic peptide from mouse PCSK9 amino acids 152-163 (VFAQSIPWNLER). This peptide is conserved in human and rat PCSK9.
<b>Host:</b>	Rabbit
<b>Cross Reactivity:</b>	(+) Human, mouse, and rat PCSK9; other species not tested
<b>Stability:</b>	≥1 year at -20°C
<b>Applications:</b>	The recommended starting dilution for western blot is <i>lot specific: lot specific</i> and 1:50 (8 µg/ml) for immunocytochemistry. Other applications were not attempted and therefore optimal working dilutions should be determined empirically.

Proprotein convertase subtilisin kexin 9 (PCSK9) is a member of the subtilisin serine protease family with an important role in lipoprotein metabolism.<sup>1</sup> Mutation in the PCSK9 gene is associated with autosomal dominant hypercholesterolemia which is characterized by an increase in low density lipoprotein (LDL) cholesterol levels.<sup>2</sup> PCSK9 overexpression in wild-type mice doubles the plasma total cholesterol, possibly through acceleration of the degradation of the LDL receptor.<sup>1,3</sup> PCSK9 mRNA is detected in various tissues such as liver, kidney, lung, spleen, jejunum, ileum, colon, and muscle with the highest expression in the liver.<sup>4</sup> Human PCSK9 precursor is 692 amino acid in length with an estimated molecular weight of 74 kDa. This proprotein is self-cleaved to form a mature protein at around 63 kDa in the golgi.<sup>5</sup>

### Laboratory Procedures

Immunofluorescent staining of cultured cells

1. Grow cells in multi-well plates until confluent.
2. Wash briefly with TBS, pH 7.4.
3. Fix the cells with 1% formaldehyde in TBS, pH 7.4, for 10 minutes.
4. Wash cells three times in TBS, containing 1% Triton-X 100 (TBST), 10 minutes each.
5. Incubate cells with 10% normal serum from the same species as the host of the secondary antibody in TBST, for 30 minutes.
6. Incubate cells with 8 µg/ml PCSK9 polyclonal antibody (recommended starting dilution; optimal dilution to be determined by end user) for one hour at room temperature.
7. Wash cells three times in TBST, 10 minutes each.
8. Incubate cells for one hour with an fluorochrome-conjugated secondary antibody at a concentration recommended by the provider.
9. Wash cells three times in TBST, 10 minutes each.
10. Examine the staining under a florescent microscope with the appropriate filter. Store the plate at 4°C in the dark for later analysis.

**WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY. NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.**

#### MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent via email to your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman warrants only to the original customer that the material will meet our specifications at the time of delivery.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have any obligation or liability, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees.

Buyer's exclusive remedy and Cayman's sole liability hereunder shall be limited to a refund of the purchase price, or at Cayman's option, the replacement, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our Warranty and Limitation of Remedy located on our website and in our catalog.

Copyright Cayman Chemical Company, 12/07/2011

### Cayman Chemical

#### Mailing address

1180 E. Ellsworth Road  
Ann Arbor, MI  
48108 USA

#### Phone

(800) 364-9897  
(734) 971-3335

#### Fax

(734) 971-3640

#### E-Mail

custserv@caymanchem.com

#### Web

www.caymanchem.com

## References

1. Maxwell, K.N., Fisher, E.A., and Breslow, J.L. Overexpression of PCSK9 accelerates the degradation of the LDLR in a post-endoplasmic reticulum compartment. *Proc. Natl. Acad. Sci. USA* **102(6)**, 2069-2074 (2005).
2. Abifadel, M., Varret, M., Rabès, J.-P., *et al.* Mutations in PCSK9 cause autosomal dominant hypercholesterolemia. *Nature Genet.* **34(2)**, 154-156 (2003).
3. Maxwell, K.N. and Breslow, J.L. Adenoviral-mediated expression of Pcsk9 in mice results in a low-density lipoprotein receptor knockout phenotype. *Proc. Natl. Acad. Sci. USA* **101(18)**, 7100-7105 (2004).
4. Seidah, N.G., Benjannet, S., Wickham, L., *et al.* The secretory proprotein convertase neural apoptosis-regulated convertase 1 (NARC-1): Liver regeneration and neuronal differentiation. *Proc. Natl. Acad. Sci. USA* **100(3)**, 928-933 (2003).
5. Maxwell, K.N. and Breslow, J.L. Proprotein convertase subtilisin kexin 9: The third locus implicated in autosomal dominant hypercholesterolemia. *Curr. Opin. Lipidol.* **16**, 167-172 (2005).

## Related Products

For a list of related products please visit: [www.caymanchem.com/catalog/10008811](http://www.caymanchem.com/catalog/10008811)

## Cayman Chemical

### Mailing address

1180 E. Ellsworth Road  
Ann Arbor, MI  
48108 USA

### Phone

(800) 364-9897  
(734) 971-3335

### Fax

(734) 971-3640

### E-Mail

[custserv@caymanchem.com](mailto:custserv@caymanchem.com)

### Web

[www.caymanchem.com](http://www.caymanchem.com)