

Product Information



PSD95 Polyclonal Antibody

Item No. 10009506

Synonym:	Postsynaptic Density Protein of 95
Supplied as:	100 µl affinity-purified antibody in 10 mM HEPES, pH 7.5, containing 150 mM NaCl, 100 µg/ml BSA, and 50% glycerol.
Host:	Rabbit
Antigen:	Peptide corresponding to amino acid residues from the N-terminal region of rat PSD95
Cross Reactivity:	(+) Rat and mouse PSD95; expected to react with bovine, human, non-human primates, and zebrafish PSD95
Stability:	≥1 year at -20°C
Application:	The recommended starting dilution for western blot is 1:1,000.

PSD95 is a very prominent component of the postsynaptic densities of synapses. It contains three PDZ domains which play key roles in its interactions with other proteins in the synapse. It has been proposed that these PDZ domains organize glutamate receptors and their associated signaling proteins and determine the size and strength of synapses.¹ Recent work suggests that interaction of the NMDAR with PSD95 *via* these PDZ domains can be regulated by phosphorylation.²

References

1. Ye, B., Liao, D.Z., Zhang, X.Q., Zhang, P.S., Dong, H.L., Haganir, R.L. GRASP-1: A neuronal RasGEF associated with the AMPA receptor/GRIP complex. *Neuron* **26**, 603-617 (2000).
2. Scannevin R.H., Haganir R.L. Postsynaptic organization and regulation of excitatory synapses. *Nat Rev Neurosci* **1**, 133-141 (2000).

Related Products

For a list of related products please visit: www.caymanchem.com/catalog/10009506

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

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, 12/14/2011