

Product Information



Dynamin (Phospho-Ser⁷⁷⁸) Polyclonal Antibody

Catalog No. 10009786

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:	Sheep
Antigen:	Phosphopeptide corresponding to amino acid residues surrounding Phospho-Ser ⁷⁷⁸ of rat dynamin
Cross Reactivity:	(+) rat dynamin; expected to react with bovine, canine, chicken, human, murine, non-human primate dynamin based on the fact that these species have 100% homology with the amino acid sequence used as antigen.
Stability:	≥1 year at -20°C
Application:	The recommended starting dilution for western blot is 1:1,000.

Dynamin is a member of a group of nerve terminal proteins called dephosphins that regulate synaptic vesicle endocytosis.¹⁻³ Cyclin dependent protein kinase 5 phosphorylates dynamin at Ser⁷⁷⁴ and Ser⁷⁷⁸.⁴ Phosphorylation of these sites on dynamin is thought to play a key role in synaptic vesicle trafficking.

References

1. Cousin, M., Tan, T.C., Robinson, P.J. Protein phosphorylation is required for endocytosis in nerve terminals: potential role for the dephosphins dynamin I and synaptojanin, but not AP180 or amphiphysin. *J. Neurochem* **76**, 105-116 (2001).
2. Graham, M.E., O'Callaghan, D.W., McMahon, H.T., *et al.* Dynamin-dependent and dynamin-independent processes contribute to the regulation of single vesicle release kinetics and quantal size. *Proc Natl Acad Sci USA* **99**, 7124-7129 (2002).
3. Tsuboi, T., Terakawa S., Scalettar, B.A., *et al.* Sweeping model of dynamin activity - Visualization of coupling between exocytosis and endocytosis under an evanescent wave microscope with green fluorescent proteins. *J. Biol. Chem.* **277**, 15957-15961 (2002).
4. Tan, T.C., Valova, V.A., Malladi, C.S., *et al.* Cdk5 is essential for synaptic vesicle endocytosis. Molecular characterization of the family of the N-methyl-D-aspartate receptor subunits. *Nat. Cell Biol.* **5**, 701-710 (2003).

Related Product

Dynamin (Phospho-Ser⁷⁷⁴) Polyclonal Antibody - Cat. No. 10009785

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, 09/08/2010

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