

PRODUCT INFORMATION



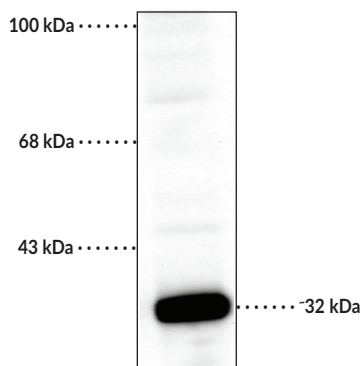
DARPP-32 Polyclonal Antibody

Item No. 29259

Overview and Properties

Contents:	This vial contains 100 µl of polyclonal antibody prepared from pooled rabbit serum via chromatography on an affinity column prepared with the N-terminal peptide used as antigen.
Synonym:	Protein Phosphatase 1 Regulatory Subunit 1B
Immunogen:	Peptide corresponding to amino acid residues from the N-terminal region of rat DARPP-32 conjugated to KLH
Molecular Weight:	~32 kDa
Species Reactivity:	(+) Mouse, rat
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥1 year
Storage Buffer:	10 mM HEPES, pH 7.5, with 150 mM sodium chloride, 100 µg/ml BSA, and 50% glycerol
Host:	Rabbit
Applications:	Immunohistochemistry (IHC) and Western blot (WB); the recommended starting dilution is 1:250 for IHC and 1:1,000 for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Image



WB of rat hippocampal lysate showing specific immunolabeling of the ~32 kDa DARPP protein. DARPP-32 Polyclonal Antibody was used as the detection antibody.

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 12/11/2020

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

Dopamine and cAMP-regulated phosphoprotein of molecular weight 32 kDa (DARPP-32) is a member of the protein phosphatase inhibitor 1 family and is encoded by the *PPP1R1B* gene in humans.¹ It is primarily expressed in medium spiny neurons of dopamine-innervated brain regions and is localized to the cytoplasm.² DARPP-32 contains multiple phosphorylation sites that allow it to function as a central regulator and integrator of signaling cascades induced by a variety of neurotransmitters, neuropeptides, and psychostimulatory agents. It also contains a protein phosphatase 1 (PP1) binding domain.³ DARPP-32 inhibits PP1 when phosphorylated at threonine 34 (Thr³⁴) by PKA and inhibits PKA when phosphorylated at Thr⁷⁵ by cyclin-dependent kinase 5 (Cdk5). *Ppp1r1b*^{-/-} mice have decreased amphetamine-induced GABA and dopamine release in the striatum and are less susceptible to catalepsy induced by raclopride (Item No. 17422) compared to wild-type mice.⁴ Increased DARPP-32 protein levels have been identified in breast, prostate, gastric, and colon cancer tumors.⁵ DARPP-32 levels are also increased in postmortem-derived dorsolateral prefrontal cortex from individuals with schizophrenia or bipolar disorder.⁶ Cayman's DARPP-32 Polyclonal Antibody can be used for immunohistochemistry (IHC) and Western blot (WB) applications. The antibody recognizes DARPP-32 at ~32 kDa from mouse and rat samples.

References

1. Meyer-Lindenberg, A., Straub, R.E., Lipska, B.K., *et al.* Genetic evidence implicating DARPP-32 in human frontostriatal structure, function, and cognition. *J. Clin. Invest.* **117**(3), 672-682 (2007).
2. Yger, M. and Girault, J.A. DARPP-32, jack of all trades... master of which? *Front. Behav. Neurosci.* **5**, 56 (2011).
3. Greengard, P., Allen, P.B., and Nairn, A.C. Beyond the dopamine receptor: The DARPP-32/protein phosphatase-1 cascade. *Neuron.* **23**(3), 435-447 (1999).
4. Fienberg, A.A., Hiroi, N., Mermelstein, P.G., *et al.* DARPP-32: Regulator of the efficacy of dopaminergic neurotransmission. *Science* **281**(5378), 838-842 (1998).
5. Belkhiri, A., Zhu, S., and El-Rifai, W. DARPP-32: From neurotransmission to cancer. *Oncotarget* **7**(14), 17631-17640 (2016).
6. Ishikawa, M., Mizukami, K., Iwakiri, M., *et al.* Immunohistochemical and immunoblot analysis of dopamine and cyclic AMP-regulated phosphoprotein, relative molecular mass. *Prog. Neuropsychopharmacol. Biol. Psychiatry* **31**(6), 1177-1181 (2007).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM