

PRODUCT INFORMATION



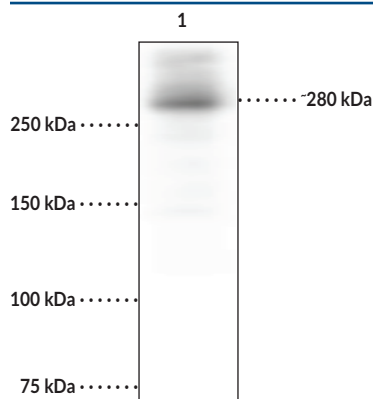
MAP2 (bovine) Monoclonal Antibody (Clone 4H5)

Item No. 29280

Overview and Properties

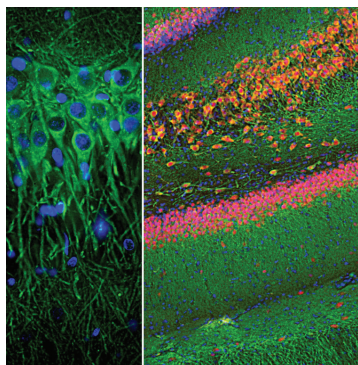
Contents:	This vial contains 100 µl of protein G-purified monoclonal antibody.
Synonyms:	MAP-2, Microtubule-associated Protein 2
Immunogen:	Purified bovine MAP2
Molecular Weight:	~280 kDa
Species Reactivity:	(+) Human, bovine, mouse, rat
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥1 year
Storage Buffer:	PBS, pH 7.2, with 5 mM sodium azide and 50% glycerol
Host:	Mouse
Applications:	Immunocytochemistry (ICC), immunohistochemistry (IHC), and Western blot (WB); the recommended starting dilution is 1:1,000. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Images



Lane 1: Mouse cortical lysate

WB of mouse cortical lysate showing specific immunolabeling of the ~280 kDa MAP2 protein.



Immunolabeling of rat hippocampus showing specific labeling of neuronal dendrites and perikarya with MAP2 (bovine) Monoclonal Antibody (Clone 4H5) in green, FOX3 in red, and nuclei in blue.

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, 12/10/2020

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

Microtubule-associated protein 2 (MAP2) belongs to the microtubule-associated protein (MAP) family of cytoskeletal filament proteins, which includes MAP1-5 and tau.¹ MAP2 is composed of an N-terminal projection domain that interacts with PKA and a C-terminal domain that contains microtubule binding repeats containing a KXGS motif that is subject to phosphorylation.² Five isoforms of MAP2 are generated through alternative splicing and are found in the cell body and dendrites of neurons, with some isoforms also found in astrocytes and oligodendrocytes of the CNS and in the testis.^{1,2} MAP2 is expressed in mature neurons, with the MAP2a and MAP2c isoforms expressed in a developmental stage-dependent manner while the MAP2b isoform is expressed throughout development and adulthood.^{1,3} MAP2 binds to and stabilizes microtubules and increases their rigidity, as well as interacts with F-actin during neurite initiation and with proteins involved in signal transduction.² Phosphorylation of the KXGS motif by MAP/microtubule affinity-regulating kinase 1 (MARK1), MARK2, or PKA decreases the affinity of MAP2 for microtubules, and the detachment of MAP2 leads to microtubule destabilization and disassembly. Decreased levels of MAP2 protein have been found in postmortem brain tissue from patients with a Lewy body variant of Alzheimer's disease.⁴ MAP2 has been found in cytoplasmic neuronal Lewy bodies colocalized with α -synuclein and ubiquitin in postmortem substantia nigra from patients with Parkinson's disease.⁵ Cayman's MAP2 (bovine) Monoclonal Antibody (Clone 4H5) can be used for immunocytochemistry (ICC), immunohistochemistry (IHC), and Western blot (WB) applications. The antibody recognizes MAP2 at approximately 280 kDa from human, bovine, mouse, and rat samples.

References

1. Kindler, S. and Garner, C.C. Four repeat MAP2 isoforms in human and rat brain. *Brain Res. Mol. Brain Res.* **26(1-2)**, 218-224 (1994).
2. Dehmelt, L. and Halpain, S. The MAP2/Tau family of microtubule-associated proteins. *Genome Biol.* **6(1)**, 204 (2004).
3. Cáceres, A., Banker, G.A., and Binder, L. Immunocytochemical localization of tubulin and microtubule-associated protein 2 during the development of hippocampal neurons in culture. *J. Neurosci.* **6(3)**, 714-722 (1986).
4. Melková, K., Zapletal, V., Narasimhan, S., *et al.* Structure and functions of microtubule associated proteins tau and MAP2c: Similarities and differences. *Biomolecules* **9(3)**, 105 (2019).
5. D'Andrea, M.R., Ilyin, S., and Plata-Salaman, C.R. Abnormal patterns of microtubule-associated protein-2 (MAP-2) immunolabeling in neuronal nuclei and Lewy bodies in Parkinson's disease substantia nigra brain tissues. *Neurosci. Lett.* **306(3)**, 137-140 (2001).

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