

PRODUCT INFORMATION



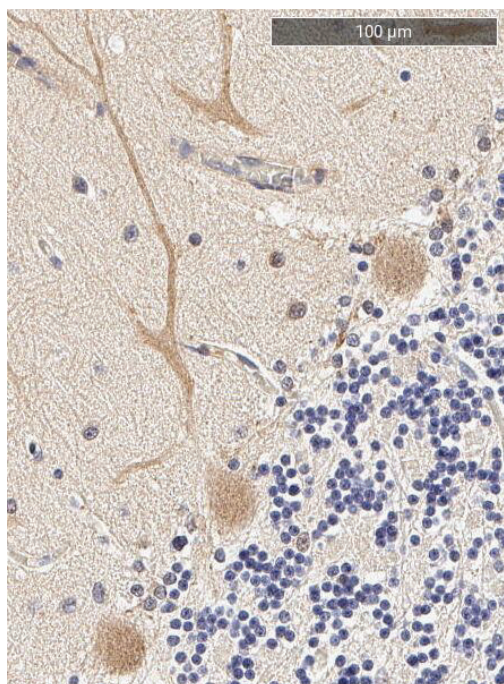
Optineurin (C-Term) Polyclonal Antibody

Item No. 100000

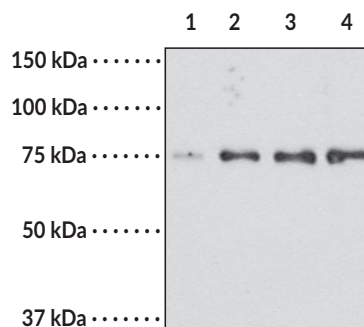
Overview and Properties

Contents:	This vial contains 500 µl of peptide affinity-purified IgG.
Synonyms:	FIP-2 (<i>14.7K interacting protein-2</i>); NRP (<i>nemo-related protein</i>)
Immunogen:	Synthetic peptide from C-terminal region of human optineurin
Species Reactivity:	(+) Human, mouse, and rat optineurin; other species not tested
Uniprot No.:	Q96CV9
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥1 year
Storage Buffer:	PBS, pH 7.2 with 50% glycerol and 0.02% sodium azide
Host:	Rabbit
Isotype:	IgG
Application:	Immunohistochemistry (IHC) and Western blot (WB); the recommended starting dilution for IHC is 1:80 and 1:200 for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Images



Immunohistochemistry analysis of formalin-fixed, paraffin-embedded (FFPE) human cerebellum tissue after heat induced antigen retrieval in pH 6.0 citrate buffer. After incubation with optineurin (C-term) polyclonal antibody (Item No. 100000) at a 1:80 dilution, slides were incubated with biotinylated secondary antibody, followed by alkaline phosphatase-streptavidin and chromogen (DAB).



Lane 1: BJ fibroblast lysates (7.5 µg)
Lane 2: HEK293 cell lysates (20 µg)
Lane 3: HEK293 cell lysates (25 µg)
Lane 4: HEK293 cell lysates (30 µg)

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, 09/14/2018

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

Optineurin, originally named FIP-2, is a 74 kDa protein implicated in signal transduction of the tumor necrosis factor (TNF) pathway as well as membrane traffic regulation and cellular morphogenesis. Optineurin protein is also involved in apoptosis, although the exact role is currently unknown. Increased apoptosis of trabecular meshwork cells is a factor in the pathogenesis of glaucoma and mutations in the optineurin gene have been correlated with the some types of the disease. Cayman Chemical is a leading provider of prostaglandin analogs, such as latanoprost and fluprostenol, currently utilized for lowering the intraocular pressure (IOP) associated with glaucoma. Our optineurin polyclonal antibody is another research tool useful for delineating the physiology of the disease.

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