

PRODUCT INFORMATION



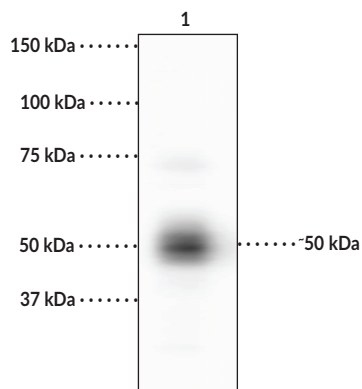
c-Fos Monoclonal Antibody

Item No. 29262

Overview and Properties

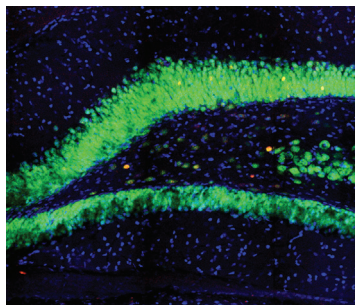
| | |
|----------------------------|---|
| Contents: | This vial contains 100 µl of protein G-purified monoclonal antibody. |
| Synonyms: | AP-1 Transcription Factor Subunit, FBJ Murine Osteosarcoma Viral Oncogene Homolog, Fos Proto-oncogene, Proto-oncogene c-Fos |
| Immunogen: | Recombinant human full-length c-Fos |
| Molecular Weight: | ~50 kDa |
| Species Reactivity: | (+) Human, bovine, mouse, rat |
| Form: | Liquid |
| Storage: | -20°C (as supplied) |
| Stability: | ≥1 year |
| Storage Buffer: | PBS 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 for ICC and IHC and 1:2,000 for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically. |

Images



Lane 1: HeLa cell lysate

WB of HeLa cell lysate showing specific immunolabeling of the ~50 kDa Fos protein. c-Fos Monoclonal Antibody was used as the detection



Immunofluorescence of a section of rat hippocampus showing neurons labeled with c-Fos Monoclonal Antibody (red), counterstained with FOX3/NeuN (green), and the nuclei stained with DAPI. When both Fos and FOX3 are expressed, the label appears orange.

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

c-Fos is a widely expressed nuclear protein and member of the Fos protein family that includes c-Fos, FosB, Fra-1, and Fra-2.¹ It is a proto-oncogene that contains a leucine zipper motif to facilitate heterodimerization with Jun family proteins, such as c-Jun, to form the transcription factor complex activator protein-1 (AP-1) that has roles in cell proliferation, differentiation, and apoptosis. Expression of *c-fos* is induced in response to growth factors, tumor promoters, cytokines, and UV radiation.^{1,2} In neurons, *c-fos* expression is induced upon depolarization and is widely used as a marker of neuronal activation.³ *c-fos* is also overexpressed in a variety of tumor cells, and transgenic overexpression of *c-fos* in mice induces formation of chondrosarcomas.^{2,4} Cayman's c-Fos Monoclonal Antibody can be used for immunocytochemistry (ICC), immunohistochemistry (IHC), and Western blot (WB) applications. The antibody recognizes c-Fos at approximately 50 kDa from human, bovine, mouse, and rat samples.

References

1. Kovács, K.J. c-Fos as a transcription factor: A stressful (re)view from a functional map. *Neurochem. Int.* **33(4)**, 287-297 (1998).
2. Tkach, V., Tulchinsky, E., Lukanidin, E., *et al.* Role of the fos family members, c-Fos, Fra-1 and Fra-2, in the regulation of cell motility. *Oncogene* **22(32)**, 5045-5054 (2003).
3. Hoffman, G.E., Smith, M.S., and Verbalis, J.G. c-Fos and related immediate early gene products as markers of activity in neuroendocrine systems. *Front. Neuroendocrinol.* **14(3)**, 173-213 (1993).
4. Jacenko, O. *c-fos* and bone loss: A proto-oncogene regulates osteoclast lineage determination. *BioEssays* **17(4)**, 277-281 (1995).

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