

PRODUCT INFORMATION



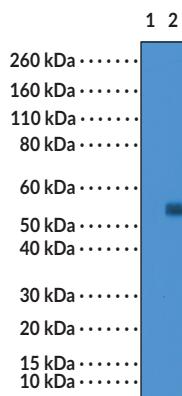
c-Fos (N-Term) Rabbit Monoclonal Antibody (Clone RM374)

Item No. 32299

Overview and Properties

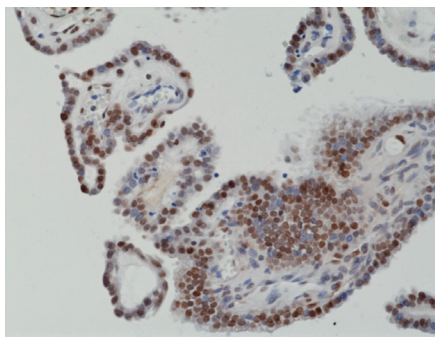
Contents:	This vial contains 100 µl of protein A-affinity purified monoclonal antibody.
Synonyms:	Cellular Oncogene Fos, Proto-oncogene c-Fos
Immunogen:	Peptide from the N-terminal region of human c-Fos
Cross Reactivity:	(+) c-Fos
Species Reactivity:	(+) Human
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥1 year
Storage Buffer:	PBS with 50% glycerol, 1% BSA, and 0.09% sodium azide
Clone:	RM374
Host:	Rabbit
Isotype:	IgG
Applications:	Immunohistochemistry (IHC) and Western blot (WB); the recommended starting dilution is 1:500-1:1,000 for IHC and 1:1,000-1:5,000 for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Images



Lane 1: HeLa cell lysates untreated
Lane 2: HeLa cell lysates treated

WB of HeLa cell lysates untreated or treated with TPA using c-Fos (N-Term) Rabbit Monoclonal Antibody (Clone RM374) at a dilution of 1:5,000.



Immunohistochemical staining of formalin-fixed and paraffin-embedded human thyroid tissue using c-Fos (N-Term) Rabbit Monoclonal Antibody (Clone RM374) at a dilution of 1:1,250.

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, 02/20/2024

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 (N-Term) Rabbit Monoclonal Antibody (Clone RM374) can be used for immunohistochemistry (IHC) and Western blot (WB) applications.

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