

PRODUCT INFORMATION



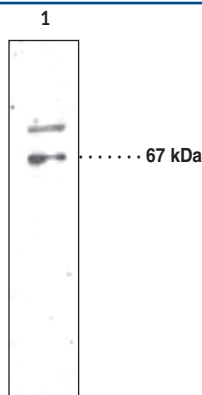
AIF Polyclonal Antibody

Item No. 160773

Overview and Properties

Contents:	This vial contains 500 μ l of peptide affinity purified polyclonal antibody.
Synonyms:	Apoptosis-Inducing Factor, Programmed Cell Death Protein 8
Immunogen:	Synthetic peptide from an internal region of human AIF
Species Reactivity:	(+) Human, rat, and mouse AIF; other species not tested
Uniprot No.:	O95831
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	\geq 3 years
Storage Buffer:	TBS, pH 7.4 with 50% glycerol, 0.5 mg/ml BSA, and 0.02% sodium azide
Host:	Rabbit
Applications:	Western blot; the recommended starting dilution is 1:200. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Images



Lane 1: A549 cell lysate (30 μ g)

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, 11/17/2023

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

Apoptosis-inducing factor (AIF) is a highly conserved mitochondrial protein with roles in redox-biochemistry and apoptosis.^{1,2} Apoptosis is a controlled process of cell death necessary for proper physiological development and maintenance. Loss of mitochondrial membrane potential results in the release of several proteins critical to the acceleration of apoptosis.³ When AIF is released from the mitochondrial intermembrane space it migrates to the nucleus to initiate chromatin condensation and DNA cleavage.⁴⁻⁶ AIF is recognized by immunoblotting at 67 kDa in most tissues and cell lines.

References

1. Lorenzo, H.K., Susin, S.A., Penninger, J., *et al.* Apoptosis inducing factor (AIF): A phylogenetically old, caspase-independent effector of cell death. *Cell Death and Differentiation* **6**, 516-524 (1999).
2. Joza, N., Susin, S.A., Daugas, E., *et al.* Essential role of the mitochondrial apoptosis-inducing factor in programmed cell death. *Nature* **410**, 549-554 (2001).
3. Green, D.R. and Reed, J.C. Mitochondria and apoptosis. *Science* **281**, 1309-1312 (1998).
4. Susin, S.A., Lorenzo, H.K., Zamzami, N., *et al.* Molecular characterization of mitochondrial apoptosis-inducing factor. *Nature* **397**, 441-446 (1999).
5. Susin, S.A., Zamzami, N., Castedo, M., *et al.* Bcl-2 inhibits the mitochondrial release of an apoptogenic protease. *J. Exp. Med.* **184**, 1331-1341 (1996).
6. Daugas, E., Susin, S.A., Zamzami, N., *et al.* Mitochondrio-nuclear translocation of AIF in apoptosis and necrosis. *FASEB J.* **14**, 729-739 (2000).

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