

PRODUCT INFORMATION



Caspase-3 Monoclonal Antibody (Clone 31A893)

Item No. 13911

Overview

Contents: This vial contains 100 µg protein G-purified IgG in 200 µl PBS containing 0.05% BSA and 0.05% sodium azide.
Synonyms: Apopain, CPP32, ICE3, Yama
Immunogen: Recombinant full-length human caspase-3
Species Reactivity: (+) Human caspase-3

Properties

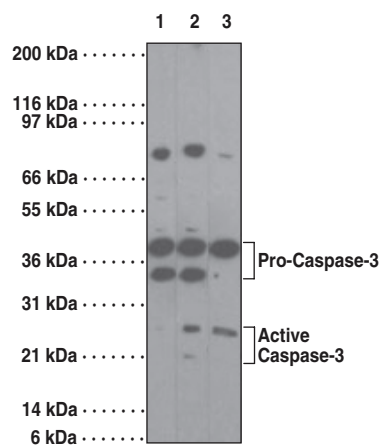
Storage: ≤6 months at 4°C; ≥6 months at -20°C
Clone: 31A893
Host: Mouse

Application

Western blot: Suggested starting concentration is 1-2 µg

Other applications were not attempted and therefore optimal working dilutions should be determined empirically.

Images



Lane 1: Jurkat cell lysate (0 hours)*
Lane 2: Jurkat cell lysate (1 hour)*
Lane 3: Jurkat cell lysate (4 hours)*

* After treatment with 2 µM staurosporine

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, 08/03/2015

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