

Caspase-2 monoclonal antibody (11B4)

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 28

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-804-356-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

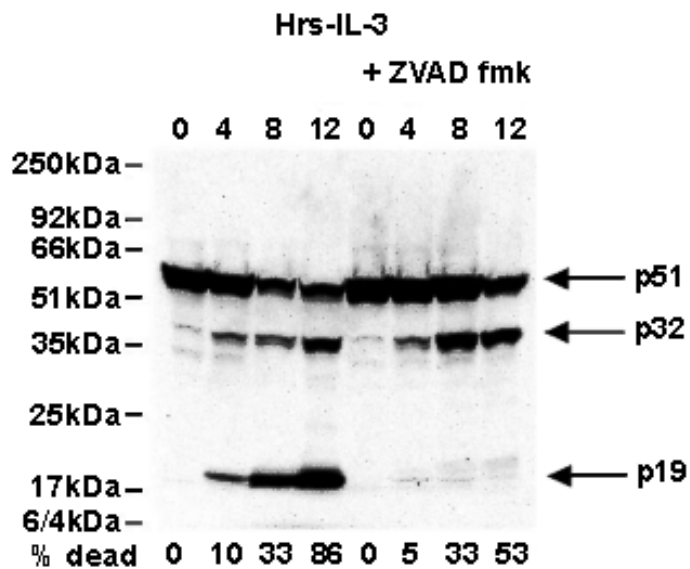
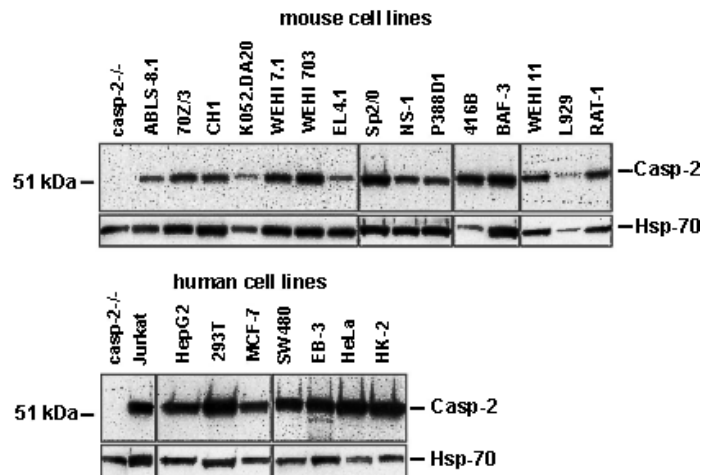


Figure: Detection of procaspase-2 (p51), an intermediate cleavage product of caspase-2 (p32) and activated caspase-2 (p19) in FDC-P1 cells (mouse IL-3 dependent promyelocytic line) during growth factor withdrawal induced apoptosis. Activation of caspase-2 is inhibited in the presence of the caspase inhibitor Z-VAD-FMK ([ALX-260-020](#) or ALX-260-138 (Ready-to-Use)).



Figures: Western blot analysis of cell lysates from human and mouse cell lines with clone 11B4. Clone 11B4 demonstrates its ability to recognize an approximately 51 kDa protein that corresponds to endogenous pro-caspase-2. Hsp-70 was used as an equal loading control, and ^{-/-} caspase-2 cells were used as a negative control.

Handling & Storage

Handling	Avoid freeze/thaw cycles.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application	IP, WB
Application Notes	Detects bands of ~51kDa (procaspase-2), ~32kDa and ~18kDa (cleavage products) by Western blot. For FC, ICC and IHC (FS) use MAb to Caspase-2 (10C6) (Prod. No. ALX-804-355).
Clone	11B4
Formulation	Liquid. In PBS containing 0.02% sodium azide.
Host	Rat
Immunogen	Recombinant human caspase-2 p19 fragment.
Isotype	IgG2a
Purity Detail	Protein G-affinity purified.
Recommendation Dilutions/Conditions	Western Blot (2µg/ml) Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.
Species Reactivity	Dog, Human, Monkey, Mouse, Rat
Specificity	Recognizes an epitope in the p19 subunit of human, mouse, rat, monkey and dog caspase-2. Detects bands of ~51kDa (procaspase-2), ~32kDa and ~18kDa (cleavage products) by Western blot.

UniProt ID

P42575

Worry-free Guarantee

This antibody is covered by our [Worry-Free Guarantee](#)

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)