

Caspase-8 (human) monoclonal antibody (12F5)

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

Citations: 80

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-804-242-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

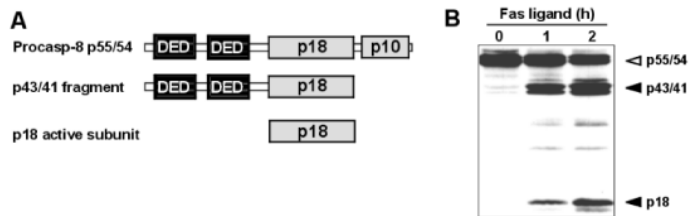


Figure:

(A) Schematic structure of caspase-8 and its cleavage products as detected by anti-caspase-8 (12F5). The antibody is directed against the p18 subunit and detects procaspase-8 (p55/54), the intermediate cleavage products of 43kD and 41kD and the p18 active subunit of caspase-8.

(B) Detection of Fas ligand-induced caspase-8 processing and activation in human Jurkat cells.

Method: Jurkat cells were treated with Fas ligand (100ng/ml). After the indicated time points cell lysates were prepared and separated on a 12.5% SDS-PAGE (2×10^6 cells/lane) under reducing conditions. Proteins were immunoblotted with anti-caspase-8 12F5 (1 μ g/ml). Following incubation with peroxidase-conjugated secondary antibodies caspase-8 processing was detected by enhanced chemoluminescent staining. The open arrowhead indicates the two isoforms of procaspase-8 (caspase-8/a and caspase-8/b), which are cleaved into the intermediate forms p43 and p41 and finally processed to the active p18 subunit (filled arrowheads).

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name FLICE

Application IHC (FS), IP, WB

Clone 12F5

Crossreactivity Does not cross-react with mouse caspase-8.

Formulation Liquid. In 0.15M PBS, pH 7.2, containing 1% BSA and 0.01% thimerosal.

Host Mouse

Immunogen Recombinant human caspase-8.

Isotype IgG2b

Purity Detail Protein G-affinity purified.

Recommendation Dilutions/Conditions Western Blot (1:100)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Species Reactivity Human

Specificity Recognizes procaspase-8 and active caspase-8. Detects procaspase-8 (p55/54), the intermediate cleavage products (p43/41) and the p18 active subunit of caspase-8 by Western blot.

UniProt ID Q14790

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)