

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

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

Citations: 78

[View Online »](#)

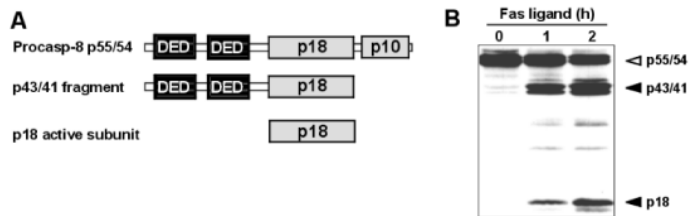
## Ordering Information

[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 ( $2 \times 10^6$  cells/lane) under reducing conditions. Proteins were immunoblotted with anti-caspase-8 12F5 (1 $\mu$ 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).

<b>Handling</b>	Avoid freeze/thaw cycles.
<b>Long Term Storage</b>	+4°C
<b>Shipping</b>	Blue Ice

## 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)