

Caspase-8 (mouse) monoclonal antibody (3B10)

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

Citations: 12

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-804-448-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

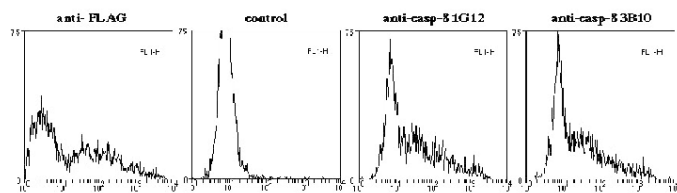


Figure 1: Flow cytometry of HEK 293T cells transfected with FLAG tagged mutated mouse caspase-8. **Method:** MAbs 1G12 and 3B10 both detect overexpressed FLAG-epitope-tagged pro-caspase-8 (active-site-mutant) by flow cytometry in permeabilized cells. As a positive control staining with a MAb to FLAG was performed. As a negative control an isotype control was employed.

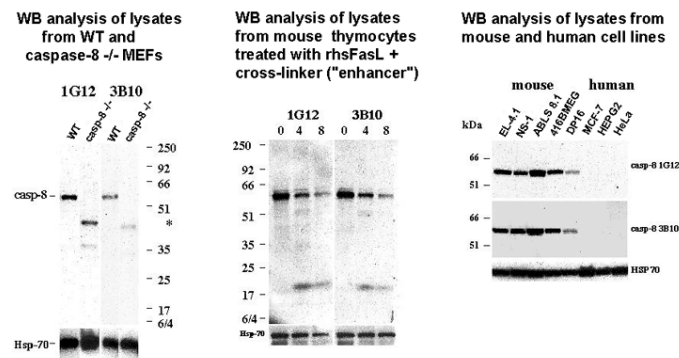


Figure 2: MAbs 1G12 and 3B10 both detect pro-caspase-8 in MEFs from WT mice, but not in MEFs from caspase-8^{-/-} mice. Several smaller bands detected in the caspase-8^{-/-} MEFs, correspond to truncated forms of caspase-8 made in the caspase-8^{-/-} mice since only exons 1 and 2 of mouse caspase-8 were deleted in these knock-out mice and not the region encoding the p18 subunit. Note: Extra bands marked by * are only seen in lysates from caspase-8^{-/-} MEFs and not in lysates from any WT cell lines or mouse WT tissue. Both MAbs to caspase-8 (mouse) do not recognize human caspase-8, whereas endogenous caspase-8 can be efficiently detected in various mouse cell lines. Upon an apoptotic stimulus e.g. by cross-linked rhsFasL both MAbs to caspase-8 (mouse) do also recognize the cleaved active p18 subunit of mouse caspase-8 in addition to the caspase-8 precursor.

Handling & Storage

Handling	For long term storage, aliquot and freeze at -20°C. 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

Alternative Name	FLICE
Application	ELISA, Flow Cytometry, ICC, WB
Application Notes	ELISA: recombinant mouse caspase-8 Flow Cytometry: overexpressed casepase-8 Excellent for Immunocytochemistry. Detects bands of ~55kDa (full-length caspase-8) and ~18kDa (apoptosis-induced cleavage fragment) by Western blot.
Clone	3B10
Crossreactivity	Does not cross-react with human caspase-8.
Formulation	Liquid. In PBS containing 0.02% sodium azide.
Host	Rat
Immunogen	Recombinant mouse caspase-8 p18 subunit.
Isotype	IgG1
Purity Detail	Protein G-affinity purified.
Species Reactivity	Mouse
UniProt ID	O89110

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

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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