

Caspase-8 (mouse) monoclonal antibody (1G12)

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

Citations: 103

[View Online »](#)

Ordering Information

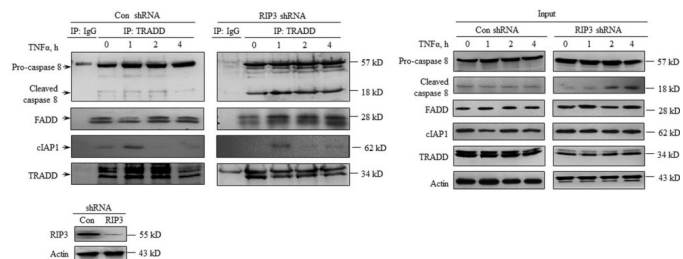
[Order Online »](#)

ALX-804-447-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

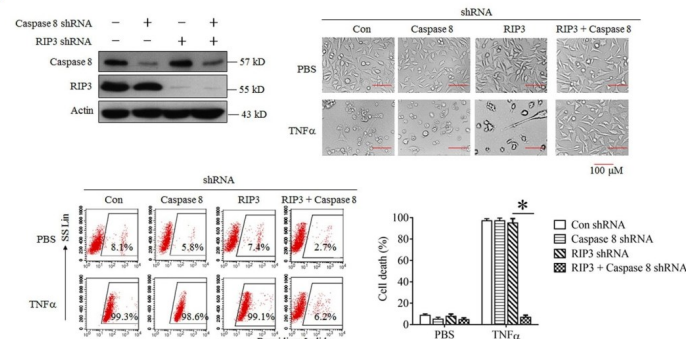
C



TRADD activates the caspase pathway by binding to and activating caspase 8. (A) TRADD mediates the TNF α -induced activation of the caspase pathway in the absence of RIP3. The negative control, TRADD knockdown, RIP3 knockdown or RIP3 and TRADD double-knockdown L929 cells were treated with TNF α for the indicated times, and the cleavage of PARP and caspase 3 was assessed by western blotting. Actin was used as a loading control. The full-length blots are presented in Supplementary Figure 4A. (B) TRADD knockdown suppresses TNF α -triggered caspase 8 activation in RIP3 knockdown L929 cells. TRADD knockdown, RIP3 knockdown or RIP3 and TRADD double-knockdown L929 cells were treated with or without TNF α for 12 h and then harvested for the measurement of caspase 8 activity. * $P < 0.01$ compared to the control shRNA group treated with TNF α . (C) RIP3 knockdown enhances the interactions between TRADD and caspase 8. RIP3 knockdown and the negative control L929 cells were treated with TNF α for the indicated times, and the cell lysates were immunoprecipitated with a TRADD antibody. Western blotting was used to detect TRADD, caspase 8, cIAP1, FADD and Actin. The full-length blots are presented in Supplementary Figure 4C.

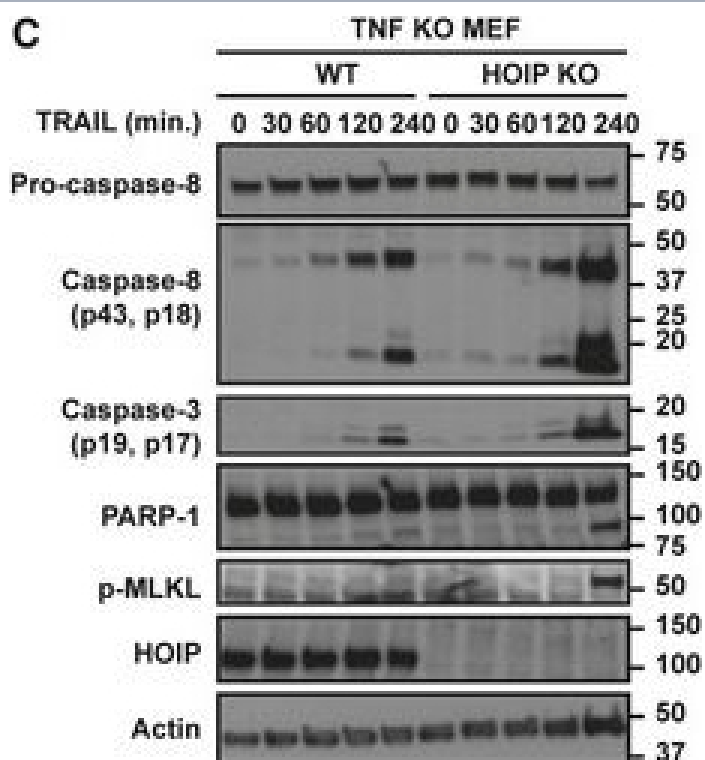
Image collected and cropped by CiteAb under a CC-BY license from the following publication: TRADD mediates the tumor necrosis factor-induced apoptosis of L929 cells in the absence of RIP3. *Sci Rep* (2017)

D



RIP3 knockdown switches TNF α -induced necroptosis to apoptosis in L929 cells. (A) Z-VAD blocks the TNF α -induced death of RIP3 knockdown L929 cells. The cells were infected with RIP3 shRNA or the control shRNA lentivirus, and western blotting was performed to determine the RIP3 knockdown efficiency. The full-length blots are presented in Supplementary Figure 1A. The cells were treated with TNF α or TNF α plus Z-VAD for 48 h, and cell death was measured by microscopy (200 \times) and flow cytometry. * $P < 0.01$. (B) RIP3 knockdown facilitates the TNF α -triggered activation of the caspase pathway. L929 cells were infected with the RIP3 shRNA or the negative control shRNA lentivirus and then treated with or without TNF α for an additional 12 h. Western blotting was performed to detect the knockdown efficiency and the cleavage of PARP and caspase 3. Actin was used as a loading control. The full-length blots are presented in Supplementary Figure 1B. (C) Caspase 8 activity was significantly increased in RIP3 knockdown L929 cells following TNF α stimulation. The RIP3 knockdown and negative control L929 cells were treated with or without TNF α for 12 h and then harvested to measure the activity of caspase 8. More than three independent experiments were performed for each group, and the relative activity of caspase 8 was calculated by normalizing the caspase 8 activity of all the groups with the activity of the negative control group. * $P < 0.01$. (D) Caspase 8 mediates the TNF α -induced death of RIP3 knockdown L929 cells. The knockdown of specific genes was mediated by infecting L929 cells with lentiviruses expressing shRNAs, and western blotting was used to evaluate the knockdown efficiency. The full-length blots are presented in Supplementary Figure 1D. The cells were treated with or without TNF α for 48 h, and cell death was measured by microscopy (200 \times) and flow cytometry. The RIP3 shRNA/DMSO and the RIP3 shRNA/TNF α FACS data presented in Figure 1D are the same as that in Figure 2A. * $P < 0.01$.

Image collected and cropped by CiteAb under a CC-BY license from the following publication: TRADD mediates the tumor necrosis factor-induced apoptosis of L929



HOIP limits TRAIL induced apoptosis and necroptosis. WT and HOIP KO TNF KO MEFs were stimulated with iz TRAIL at the indicated concentrations for 24 h (n = 5; mean ± SEM). WT and HOIP KO TNF KO MEFs, pre treated with zVAD and Nec 1s as indicated for 1 h, were stimulated with iz TRAIL for 24 h (1 µg/ml) (n = 4; mean ± SEM). WT and HOIP KO TNF KO MEFs were stimulated with iz TRAIL (1 µg/ml) for the indicated times. Lysates were analysed by Western blot. Control (CTRL) and HOIP KO K562 cells were stimulated with iz TRAIL at the indicated concentrations for 24 h (n = 4; mean ± SEM). Control and HOIP KO K562 cells, pre treated with zVAD as indicated for 1 h, were stimulated with iz TRAIL (1 µg/ml) for 24 h (n = 4; mean ± SEM). Control and HOIP deficient K562 cells were stimulated with iz TRAIL (100 ng/ml) for the indicated times. Lysates were analysed by Western blot. Black arrowhead indicates the cleaved form of HOIP. Data information: Cell death was determined after 24 h of stimulation by flow cytometry after propidium iodide (PI) labelling. *P < 0.05, **P < 0.01, ***P < 0.001; statistics were performed using t test. See also Appendix Fig S1.

Image collected and cropped by CiteAb under a CC-BY license from the following publication: The linear ubiquitin chain assembly complex regulates TRAIL-induced gene activation and cell death. *EMBO J* (2017)

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 Western blot. Detects bands of ~55kDa (full-length caspase-8) and ~18kDa (apoptosis-induced cleavage fragment) by Western blot.
Clone	1G12
Crossreactivity	Does not cross-react with human caspase-8.
Formulation	Liquid. In PBS containing 0.02% sodium azide.
Host	Rat
Immunogen	Recombinant mouse p20 caspase-8 subunit.
Isotype	IgG1
Purity Detail	Protein G-affinity purified.
Species Reactivity	Mouse
Specificity	Recognizes the p18 subunit of caspase-8.

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](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)