

FasL monoclonal antibody (3C82)

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

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-804-2730	100µg
--------------	-------

Manuals, SDS & CofA

[View Online »](#)

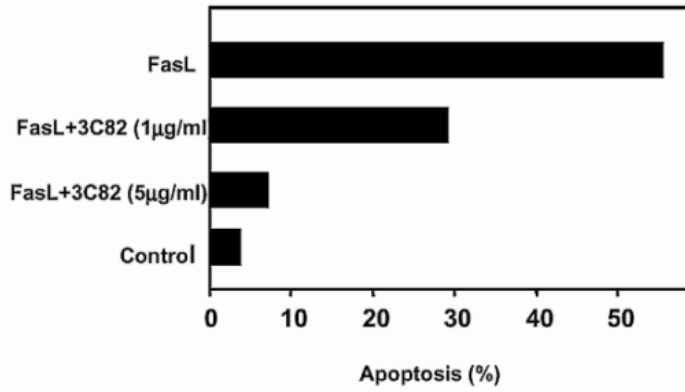
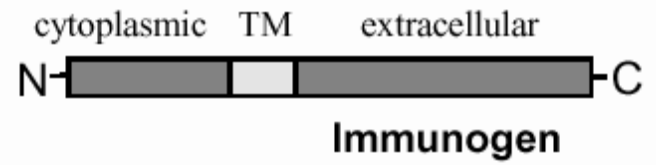


Figure: Inhibition of mouse FasL (supernatant of FasL-transfected Neuro2A cells) by 3C82. **Method:** Supernatants of Neuro2A cells containing murine FasL (1:2 dilution) were incubated with Fas expressing L929-APO-1 target cells. After 14h apoptosis was assessed by MTT staining.



Schematic structure of recombinant human FasL (extracellular domain).

Handling & Storage

Handling Do not freeze.

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name CD95L, APO-1L, CD178, TNFSF 6

Application Flow Cytometry, FUNC

Application Notes Functional Application: neutralizing

Clone 3C82

Formulation Liquid. In PBS, pH 7.2, containing 1% BSA. Contains no preservatives.

Host Mouse

Immunogen Recombinant human FasL (APO-1L; CD95L; CD178) (extracellular domain).

Isotype IgG1

Purity Detail Epitope-affinity purified.

Recommendation Dilutions/Conditions Suggested starting dilution is >1:100. Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

Species Reactivity Human, Mouse

UniProt ID P48023

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)