

Fas (human) monoclonal antibody (APO-1-1)

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

Citations: 7

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-805-022-C100	Purified - 100µg
------------------	------------------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.
Long Term Storage	+4°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	CD95, APO-1, TNFRSF 6
Application	Flow Cytometry, FUNC, IHC (FS)
Application Notes	<p>APO-1-1 (IgG1) is an isotype-switch variant of the original clone APO-1-3 (ALX-805-020). Since undesired side reactions in immunohistochemistry are minimal for IgG1 antibodies, APO-1-1 is especially useful for immunohistology on cryostat sections and on cytospin preparations.</p> <p>Not suitable for stainings on paraffin sections.</p> <p>Functional Application: Induces apoptosis upon cross-linking.</p>
Clone	APO-1-1
Formulation	Liquid. In PBS, pH 7.2-7.4.
Host	Mouse
Immunogen	Human Fas (CD95; APO-1).
Isotype	IgG1
Recommendation Dilutions/Conditions	Flow Cytometry (1-2µg/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human
UniProt ID	P25445

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)