

Fas (human) monoclonal antibody (blocking) (SM1/23)

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

Citations: 4

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-805-0100	Purified - 100µg
--------------	------------------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	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	CD95, APO-1, TNFRSF 6
Application	Flow Cytometry, FUNC
Application Notes	Functional Application: blocks induction of apoptosis induced by clone SM1/1 (Prod. No. ALX-805-008) or other commercial anti-Fas antibodies.
Clone	SM1/23
Formulation	Liquid. In PBS, pH 7.2-7.4.
Host	Mouse
Immunogen	Human Fas (CD95; APO-1).
Isotype	IgG2b
Purity Detail	Protein A affinity purified.
Species Reactivity	Human
UniProt ID	P25445

Worry-free Guarantee

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



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

Last modified: May 29, 2024