

# 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-805-010-C100	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 <a href="#">Worry-Free Guarantee</a> .



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)