

# Fas (human) monoclonal antibody (2R2)

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

Citations: 4

[View Online »](#)

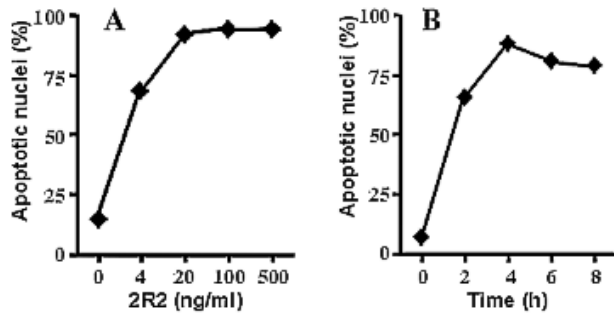
## Ordering Information

[Order Online »](#)

ALX-805-038-C100	100µg
------------------	-------

Manuals, SDS & CofA

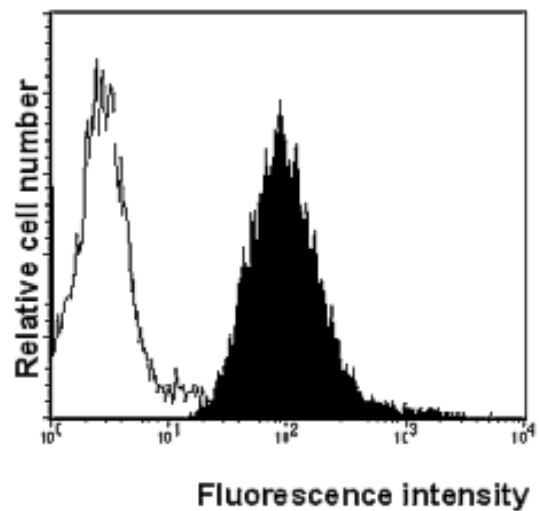
[View Online »](#)



**Figure:** Induction of apoptosis in anti-Fas 2R2 stimulated Jurkat T-cells.

**Method:** (A) Dose response: Jurkat cells ( $1 \times 10^5$ ) were treated with agonistic anti-Fas 2R2 antibodies at the indicated concentrations for 24 h. Induction of apoptosis was assessed by PI staining of hypodiploid nuclei and subsequent flow cytometry.

(B) Time course of apoptosis: Jurkat cells ( $1 \times 10^5$ ) were treated with 1  $\mu$ g/ml anti-Fas 2R2. After the indicated time points apoptosis was determined by FACS measurement of hypodiploid DNA.



**Figure:** Detection of Fas surface expression by FACS analysis.

**Method:** Normal L cells (open histogram) and cells stably overexpressing human Fas (filled histogram) were incubated with anti-Fas 2R2 and analyzed by indirect immunofluorescence using PE-labeled secondary antibodies. No immunostaining was obtained with an isotype matched IgG control.

## 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	<p>Flow Cytometry</p> <p>Functional Application (induces apoptosis – no cross-linking required).</p> <p>– Can be used in DISC (death-inducing signalling complex assays): Co-immunoprecipitates FADD and caspase-8 after receptor activation.</p> <p>Recommended working dilutions are 1:200-1:1000.</p>
Clone	2R2
Formulation	Liquid. In PBS containing 1% BSA. Contains no preservatives.
Host	Mouse
Immunogen	Mouse L cells stably overexpressing human Fas (CD95; APO-1).
Isotype	IgG3
Purity Detail	Protein A-affinity purified.

**Recommendation Dilutions/Conditions**

Flow Cytometry, Functional Application (induces apoptosis – no cross-linking required). , – Can be used in DISC (death-inducing signalling complex assays): Co-immunoprecipitates FADD and caspase-8 after receptor activation., Recommended working dilutions are 1:200-1:1,000.

**Source**

Purified from concentrated hybridoma tissue culture supernatant.

**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)