

# Fas (human) monoclonal antibody (4C3)

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

Citations: 1

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

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

**Regulatory Status** RUO - Research Use Only

## Product Details

|  |   |
|--|---|
| <b>Alternative Name</b>                    | CD95, APO-1, TNFRSF 6   |
| <b>Application</b>                         | ELISA, Flow Cytometry, WB   |
| <b>Application Notes</b>                   | Detects a band of approximately 45 kDa by Western Blot (predicted molecular weight: 37 kDa).  |
| <b>Clone</b>                               | 4C3   |
| <b>Formulation</b>                         | Liquid. In PBS. Contains no preservatives.  |
| <b>Host</b>                                | Mouse   |
| <b>Immunogen</b>                           | Human recombinant Fas (CD95; APO-1).  |
| <b>Isotype</b>                             | IgG1κ   |
| <b>Purity Detail</b>                       | Protein G affinity purified   |
| <b>Recommendation Dilutions/Conditions</b> | Flow Cytometry (1:50)Western Blot (1:1,000, ECL)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application. |
| <b>Source</b>                              | Purified from mouse ascites.  |
| <b>Species Reactivity</b>                  | Human   |
| <b>UniProt ID</b>                          | P25445  |

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)