

# Fluorescein Recombinant monoclonal antibody (4-4-20 enhanced) (Mouse Fab fragment

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

Manuals, SDS & CoFA

[View Online »](#)

## Handling & Storage

<b>Use/Stability</b>	Stable at 4°C for up to 3 months. For longer storage, aliquot and store at -20°C.
<b>Handling</b>	Avoid freeze/thaw cycles.
<b>Short Term Storage</b>	+4°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>	Fluorescyl
<b>Application</b>	Flow Cytometry, WB
<b>Application Notes</b>	Negative control
<b>Clone</b>	4-4-20 (enhanced)
<b>Formulation</b>	Liquid. In PBS containing 0.02% Proclin 300.
<b>Host</b>	Mouse
<b>Immunogen</b>	Fluorescein I-bovine serum albumin (FII-BSA).
<b>Isotype</b>	Fab fragment $\kappa$
<b>Purity Detail</b>	Immobilized metal affinity chromatography purified.
<b>Technical Info / Product Notes</b>	This is a recombinant chimeric Fab fragment created as part of a panel offering antibodies of the same specificity in different formats (species, isotype, subtype and modified versions) for use as isotype controls.
<b>Worry-free Guarantee</b>	This antibody is covered by our <a href="#">Worry-Free Guarantee</a>

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)