

# Fluorescein Recombinant monoclonal antibody (4-4-20 enhanced) (Hamster IgGκ)

This antibody is covered by our [Money-Back Guarantee](#).

## Ordering Information

[Order Online »](#)

ENZ-ABS495-0200	200µg
-----------------	-------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Use/Stability	Stable at 4°C for up to 3 months. For longer storage, aliquot and store at -20°C.
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	Fluorescyl
Application	Flow Cytometry, WB
Application Notes	Negative control
Clone	4-4-20 (enhanced)
Formulation	Liquid. In PBS containing 0.02% Proclin 300.
Host	Hamster (Armenian)
Immunogen	Fluorescein I-bovine serum albumin (FI-BSA).
Isotype	IgGk
Purity Detail	Protein A affinity purified.
Technical Info / Product Notes	This is a chimeric antibody 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.

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



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)