

# CA19-9 Recombinant monoclonal antibody (CA19.9/1390R)

CA19-9, a carbohydrate epitope expressed on a high MW (>400kDa) mucin glycoprotein, is a sialyl Lewis a structure which is synthesized from type 1 blood group precursor chains and is present in individuals expressing the Lewis a and/or Lewis b blood group antigens.

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

## Ordering Information

[Order Online »](#)

ENZ-ABS356-0100	100µg
-----------------	-------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

Application Flow Cytometry, IF, IHC (PS)

Clone CA19.9/1390R

Formulation Liquid. In 1X PBS containing 0.1 mg/ml BSA (US sourced) and 0.05% sodium azide.

Host Rabbit

Immunogen Purified human CA19-9.

Isotype IgGk

Purity Detail Protein A affinity purified.

**Recommendation** Flow Cytometry (2-4 µg/10<sup>6</sup> cells in 0.1 mL)Immunofluorescence (5-10 µg/mL)Immunohistochemistry (formalin-fixed paraffin sections) (5-10 µg/mL for 30 min at RT)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

**Dilutions/Conditions**

Species Reactivity Human

**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)