

# CEACAM1

## recombinant monoclonal antibody (103)

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

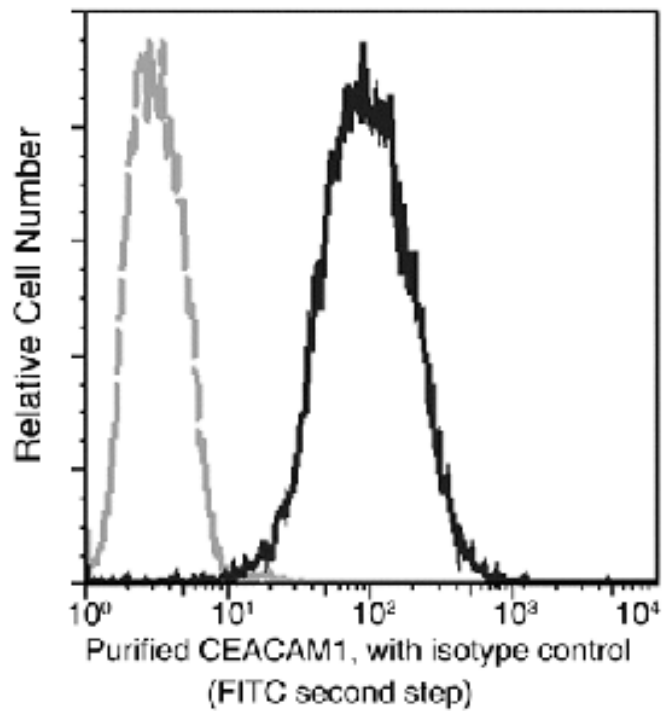
### Ordering Information

[Order Online »](#)

ENZ-ABS962-0100	100µl
-----------------	-------

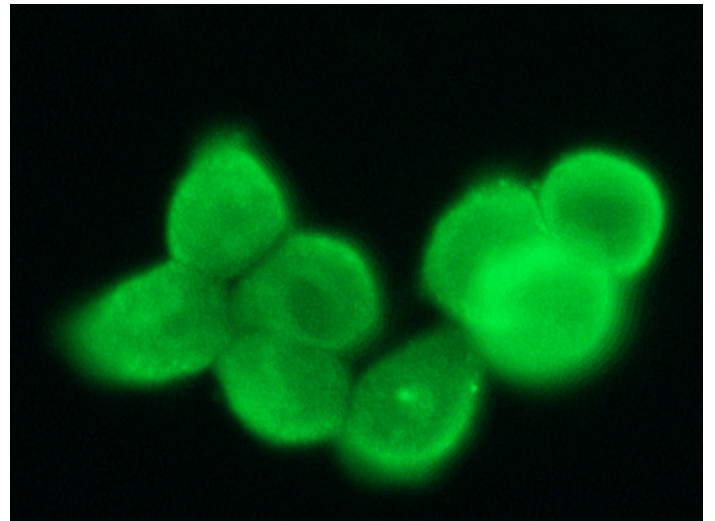
### Manuals, SDS & CofA

[View Online »](#)



© Sino Biological Inc. All rights reserved

Flow cytometric analysis of Human CEACAM1(CD66a) expression in HT29 cells. The cells were harvested and stained with purified Rabbit anti-human CEACAM1, then a FITC-conjugated second step antibody.



Immunofluorescence staining of CEACAM1 in HT29 cells. Cells were fixed with 4% PFA, permeabilized with 0.3% Triton X-100 in PBS, blocked with 10% serum, and incubated with rabbit anti-human CEACAM1 monoclonal antibody (1:60) at 4 °C overnight. Then cells were stained with the Alexa Fluor® 488-conjugated Goat Anti-rabbit IgG secondary antibody (green).

## Handling & Storage

**Handling** Protected from light. Avoid freeze-thaw cycles.

**Long Term Storage** -20°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** CD66a, Biliary glycoprotein 1, BGP

**Application** Flow Cytometry, ICC, IF

**Clone** 103

**Formulation** Liquid. 0.2 µm filtered solution in PBS. Preservative free.

**Host** Rabbit

**Immunogen** Recombinant human CEACAM1 / CD66a protein.

**Isotype** IgG

**Purity Detail** Protein A purified.

**Recommendation Dilutions/Conditions** Immunocytochemistry (1:20-1:100)Immunofluorescence (1:20-1:100)Flow Cytometry (1:25-1:100)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

**Species Reactivity** Human

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

.



Last modified: September 17, 2024

ENZO LIFE SCIENCES

Phone: 800.942.0430  
info-  
[usa@enzolifesciences.com](mailto:usa@enzolifesciences.com)

European Sales Office

ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
info-  
[eu@enzolifesciences.com](mailto:eu@enzolifesciences.com)

Belgium, The Netherlands

& Luxembourg  
Phone: +32 3 466 0420  
info-  
[be@enzolifesciences.com](mailto:be@enzolifesciences.com)

France

Phone: +33 472 440 655  
info-  
[fr@enzolifesciences.com](mailto:fr@enzolifesciences.com)

Germany

Phone: +49 7621 5500 526  
info-  
[de@enzolifesciences.com](mailto:de@enzolifesciences.com)

UK & Ireland

Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
info-  
[uk@enzolifesciences.com](mailto:uk@enzolifesciences.com)