

CEACAM1

recombinant monoclonal antibody (103)

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

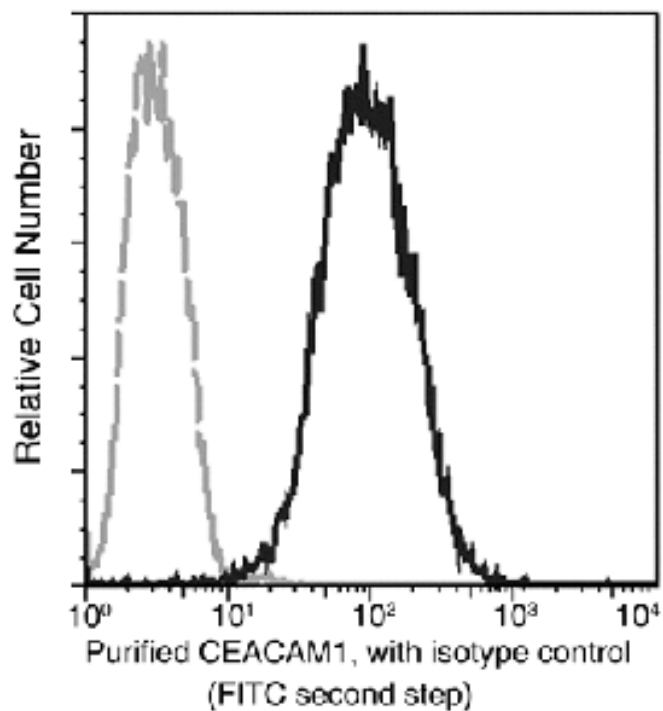
Ordering Information

[Order Online »](#)

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

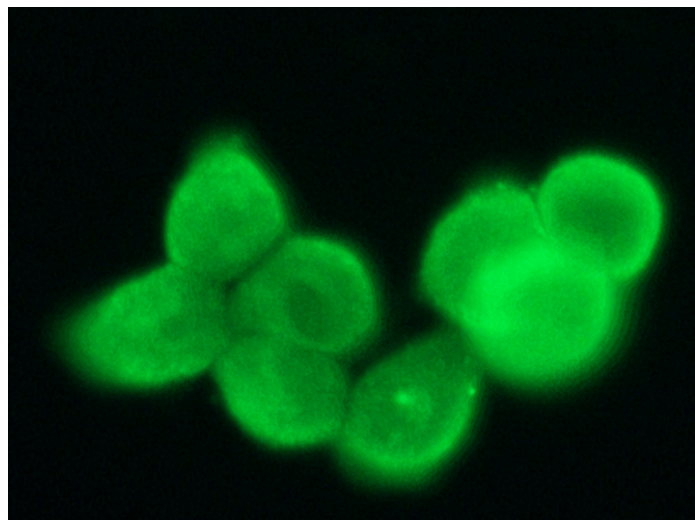
Manuals, SDS & CofA

[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 Immunocytochemistry (1:20-1:100)Immunofluorescence (1:20-1:100)Flow Cytometry
Dilutions/Conditions (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](#).



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)