

CEACAM1

recombinant monoclonal antibody (117)

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

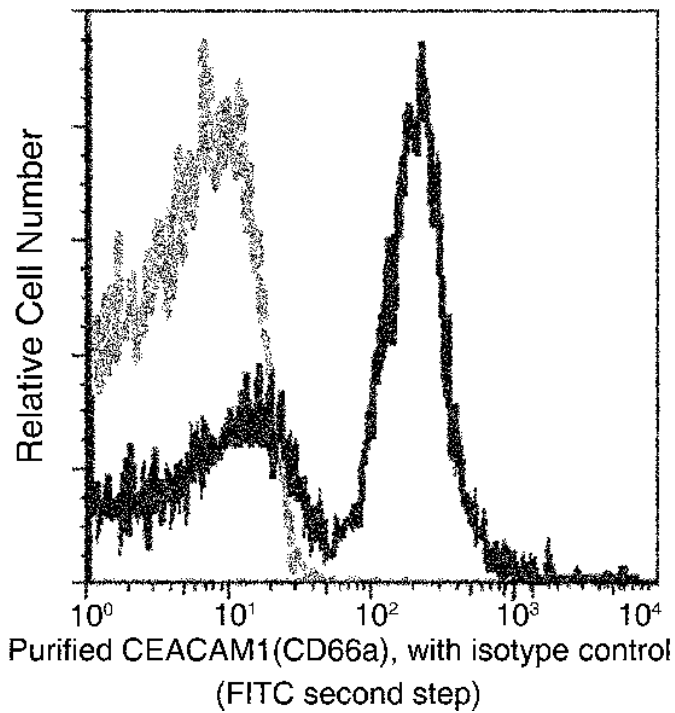
Ordering Information

[Order Online »](#)

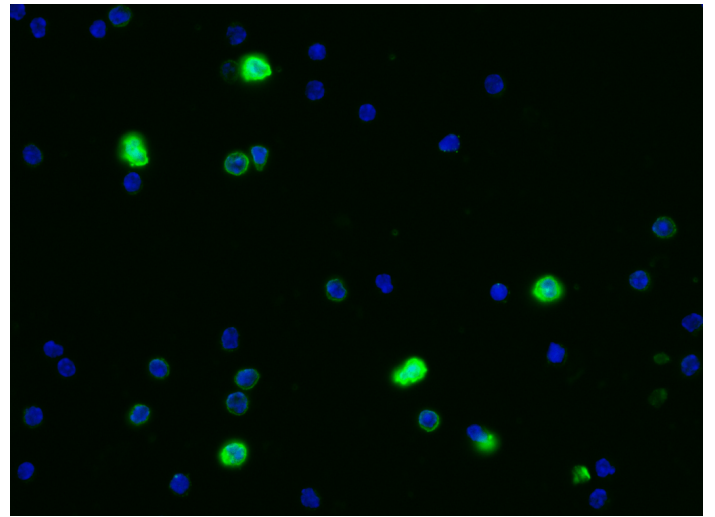
ENZ-ABS1019-0100	100µl
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Flow cytometric analysis of Mouse CEACAM1(CD66a) expression on BABL/c splenocytes. Cells were stained with purified anti-Mouse CEACAM1(CD66a), then a FITC-conjugated second step antibody. The fluorescence histograms were derived from gated events with the forward and side light-scatter characteristics of intact cells.



Immunofluorescence staining of mouse CEACAM1 in mouse splenocytes. Cells were fixed with 4% PFA, blocked with 10% serum, and incubated with rabbit anti-mouse CEACAM1 monoclonal antibody (1:60) at 4 overnight. Then cells were stained with the Alexa Fluor® 488-conjugated Goat Anti-rabbit IgG secondary antibody (green) and counterstained with DAPI (blue).

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	117
Formulation	Liquid. 0.2 µm filtered solution in PBS. Preservative free.
Host	Rabbit
Immunogen	Recombinant mouse 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	Mouse
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)