

CEACAM6

monoclonal antibody

(5E10C7)

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

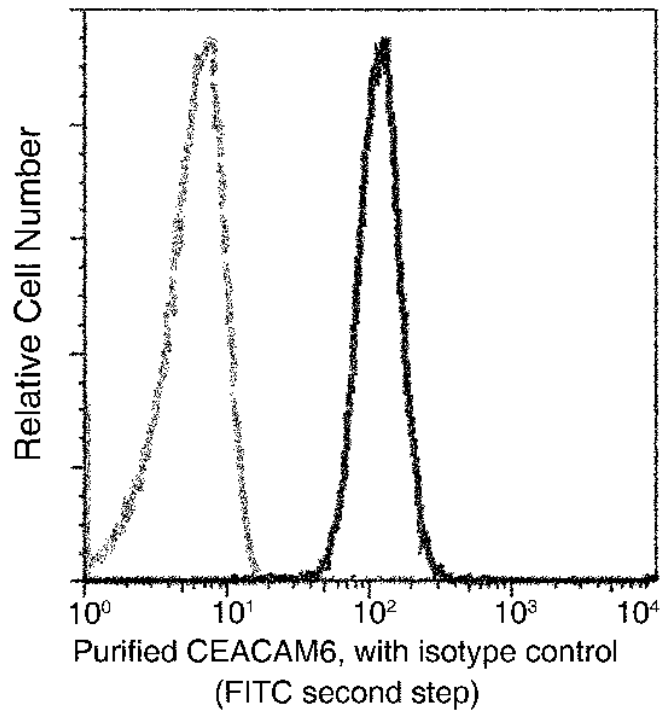
Ordering Information

[Order Online »](#)

ENZ-ABS963-0100	100µl
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Flow cytometric analysis of Human CEACAM6 expression on human whole blood granulocytes. Cells were stained with purified anti-Human CEACAM6, then a FITC-conjugated second step antibody. The histogram were derived from gated events with the forward and side light-scatter characteristics of viable granulocytes.

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 CD66c, NCA

Application ELISA, Flow Cytometry

Clone 5E10C7

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

Host Mouse

Immunogen Recombinant human CD66c protein.

Isotype IgG1

Purity Detail Protein A purified.

Recommendation Dilutions/Conditions ELISA (1:1'000-1:2'000)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,
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-
fr@enzolifesciences.com](mailto:info-fr@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)