

CEACAM6 (human) monoclonal antibody (9A6)

The normal cross-reacting antigen (NCA) or CEACAM6 belongs to the carcinoembryonic antigen (CEA) gene family. CEACAM6 is a GPI-anchored membrane protein with a single N domain, which shows structural homology to the immunoglobulin variable domains, followed by two immunoglobulin constant-like domains.

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

Citations: 15

[View Online »](#)

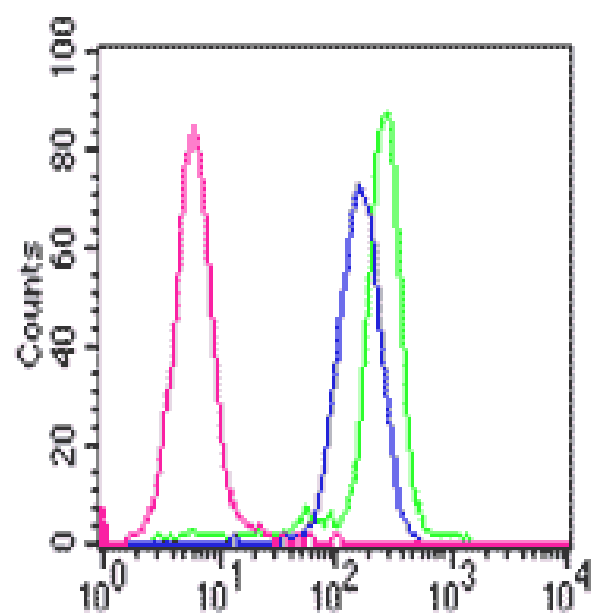
Ordering Information

[Order Online »](#)

ALX-805-083-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Handling	For maximum product recovery after thawing, centrifuge the vial before opening the cap. Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	CD66c, NCA
Application	ELISA, Flow Cytometry, IHC (FS), WB
Clone	9A6
Crossreactivity	Does not cross-react with human CEACAM1, 3, 4, 5, 7 or 8.
Formulation	Liquid. In PBS containing no preservatives.
Host	Mouse
Immunogen	Tumor cell lines expressing human CEACAM6.
Isotype	IgG1
Purity Detail	Protein G-affinity purified.
Quality Control	Routinely tested by flow cytometry on BOSC cells transiently transfected with a NCA expression vector.
Species Reactivity	Human
UniProt ID	P40199
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)