

# 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 »](#)

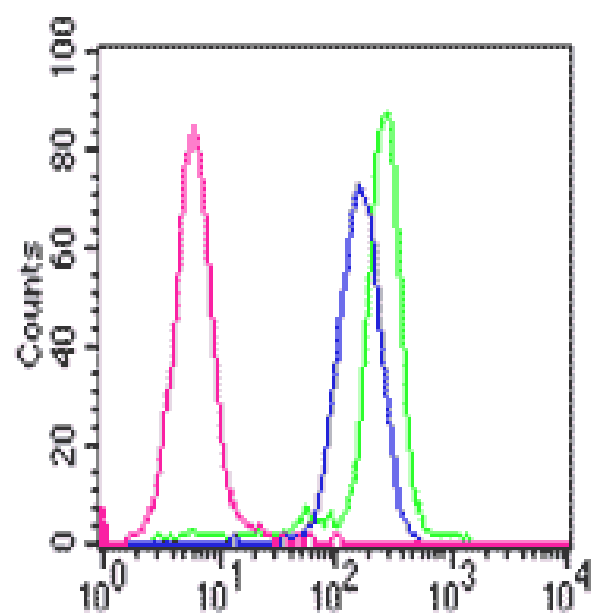
## Ordering Information

[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, pH 7.2. Contains 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



ENZO LIFE SCIENCES,  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
Phone: +32 3 466 0420  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 1 402 44 655  
[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)

This antibody is covered by our **Worry Free Guarantee**