

CEACAM1 (human) monoclonal antibody (GM8G5)

CEACAM1 (CD66a, BGP) belongs to the carcinoembryonic antigen (CEA) gene family. It is expressed on cells of epithelial and myeloid origin. Like all members of the CEACAM family, it consists of a single N domain, with structural homology to the immunoglobulin variable domains, followed by two immunoglobulin constant-like A (A1, A2) and one B domain. While N, A1 and B domains can also be found in other CEA-family members, the A2 domain of CEACAM1 differs from those found in other CEACAMs.

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

Citations: 2

[View Online »](#)

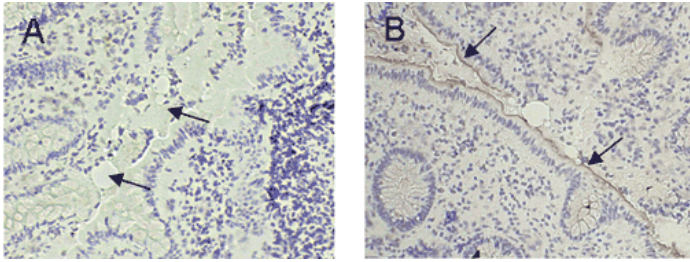
Ordering Information

[Order Online »](#)

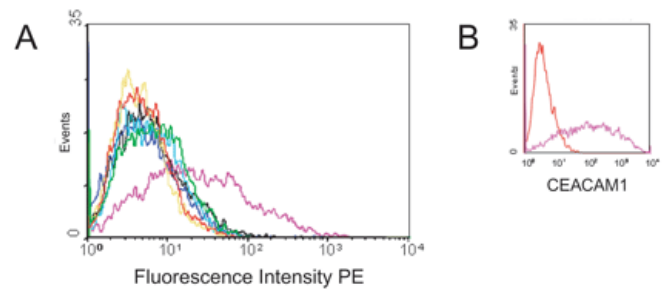
ALX-805-076-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Immunohistological staining of colon tissue with MAb to CEACAM1 (GM8G5) (Prod. No. ALX-805-076). Cryosections of colon tissue were stained with 10 $\mu\text{g}/\text{ml}$ GM8G5 diluted in PBS containing 2.5% horse serum (B). PBS 2.5% horse serum served as negative control (A). Arrows indicate representative sections of analysed tissues. Detection of GM8G5 occurred with a biotinylated anti-mouse-IgG secondary antibody and a streptavidin-peroxidase conjugate. Diaminobenzidine was used as substrate. Nuclei were stained with hematoxylin.



Specificity testing of GM8G5. BOSC cells were transiently transfected with an expression vector containing the cDNA of CEACAM1. Expression of the constructs was tested with MAb to CEACAM1 (GM8G5) (Prod. No. ALX-805-076) (B, purple curves); an irrelevant monoclonal antibody served as negative control (red curve). For specificity testing (A), GM8G5 hybridoma supernatant was tested on all CEACAM transfectants. A positive signal was obtained only with CEACAM1 transfected cells (purple curve).

Handling & Storage

Handling	For maximum product recovery after thawing, centrifuge the vial before opening the cap. Avoid freeze/thaw cycles.
Short Term Storage	+4°C
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, IHC (FS)
Clone	GM8G5
Crossreactivity	Does not cross-react with human CEACAM3, 4, 5, 6 or 8.
Formulation	Liquid. In PBS, pH 7.2. Contains no preservatives.
Host	Mouse
Immunogen	Vector containing the cDNA of human CEACAM1 A2 domain.
Isotype	IgG1
Purity Detail	Protein G-affinity purified.
Quality Control	Routinely tested on BOSC cells transiently transfected with a CEACAM1 expression vector.
Species Reactivity	Human
UniProt ID	P13688
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee

Last modified: May 29, 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-
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)