

CD66e (human) monoclonal antibody (26/3/13)

CD66e (CEACAM5; CEA) belongs to the carcinoembryonic antigen (CEA) gene family. It encodes a glycosyl phosphatidyl inositol (GPI)-linked glycoprotein with a M_r of 180'000-200'000 which is most strongly expressed on epithelial cells of the fetal and adult colon and to a minor extent on epithelial cells of the stomach and sweat glands, squamous epithelial cells of the tongue, esophagus and cervix. CD66e is used as a tumor marker for early detection of recurrent disease due to its expression in adenocarcinomas of the colon, lung, breast, stomach and pancreas and in mucinous ovarian carcinomas. Like all members of the CEACAM family, it consists of a single N domain, with structural homology to the immunoglobulin variable domains, followed by six immunoglobulin constant-like A (A1, A2, A3) and B domains (B1, B2, B3).

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

Citations: 6

[View Online »](#)

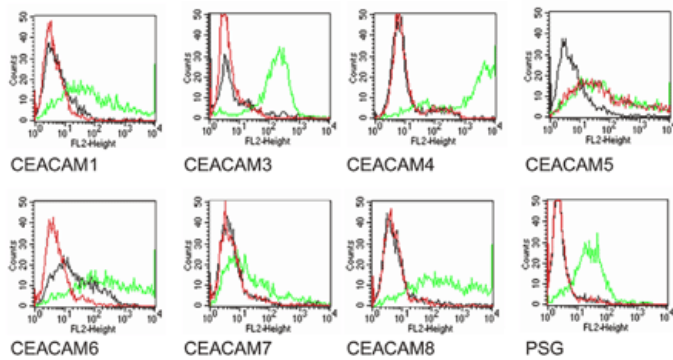
Ordering Information

[Order Online »](#)

ALX-805-077-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Specificity testing of 26/3/13 (Prod. No. ALX-805-077). Members of the CEA family were expressed on BOSC cells after transient transfection with expression vectors containing either the cDNA of CEACAM1, 5, 6, 7 or 8. Recognition of CEACAM3, 4 and of a recombinant transmembrane-anchor-ed PSG1 fusion protein was tested on stably transfected HeLa (CEACAM3, PSG) and CHO cells (CEACAM4), respectively. Expression of the constructs was confirmed with monoclonal antibodies known to recognize the corresponding proteins (CEACAM1, 3, 4, 5 and 6: D14HD11 (Prod. No. ALX-805-080); CEACAM7: CAC2; CEACAM8: TET2; PSG: BAP1; green curves). An irrelevant monoclonal antibody served as a negative control (black curves). For specificity testing, protein G purified 26/3/13 was tested on all CEACAM transfectants. A positive signal was only obtained with CEACAM5 expressing cells (red curves).

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	CEACAM5, CEA, Carcinoembryonic antigen
Application	ELISA, Flow Cytometry, IHC (FS), WB
Clone	26/3/13
Crossreactivity	Does not cross-react with human CEACAM1, 3, 4, 6, 7, 8 or PSG.
Formulation	Liquid. In PBS containing no preservatives.
Host	Mouse
Immunogen	CD66e (CEACAM5; CEA) partially purified from a perchloric acid extract from liver metastases of colonic tumors.
Isotype	IgG1
Purity Detail	Protein G-affinity purified.
Quality Control	Routinely tested by flow cytometry on BOSC cells transiently transfected with a CD66e expression vector.
Species Reactivity	Human
UniProt ID	P06731

Worry-free Guarantee

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

Last modified: December 18, 2025



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)