

CEACAM19 (human) monoclonal antibody (HY-8H10)

CEA-related cell adhesion molecule 19 (CEACAM19) belongs to the carcinoembryonic antigen (CEA) gene family (1). It encodes a putative glycoprotein which is membrane-bound via a transmembrane domain. Like all members of the CEACAM family, the CEACAM19 protein is a single-pass type I membrane protein and consists of a single extracellular N domain. Two isoforms of CEA-CAM19 exist as a result of alternative splicing events. CEACAM19 expression is ubiquitous with highest expression in prostate, uterus, fetal brain, mammary gland, adrenal gland, skeletal muscle, small intestine, and kidney. Additionally, CEACAM19 is highly expressed in breast and prostate cancer cell lines.

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

Citations: 1

[View Online »](#)

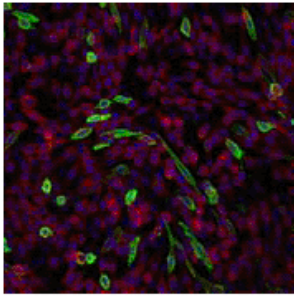
Ordering Information

[Order Online »](#)

ALX-805-0890	100µg
--------------	-------

Manuals, SDS & CofA

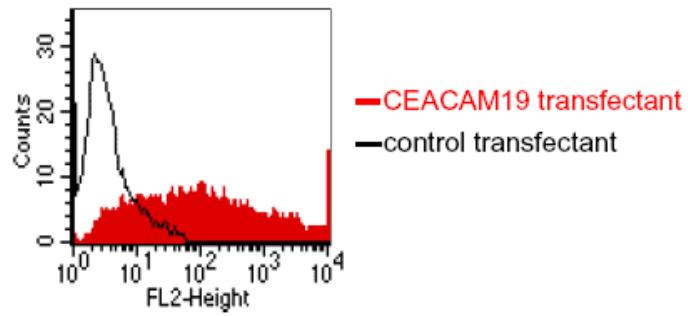
[View Online »](#)



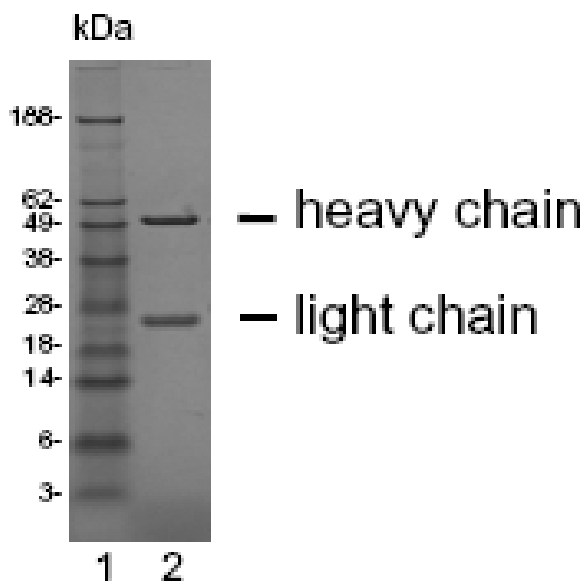
— anti-CEACAM19 (HY-8H10)
— Actin filaments
— Nuclei

Specificity testing by spectral confocal microscopy.

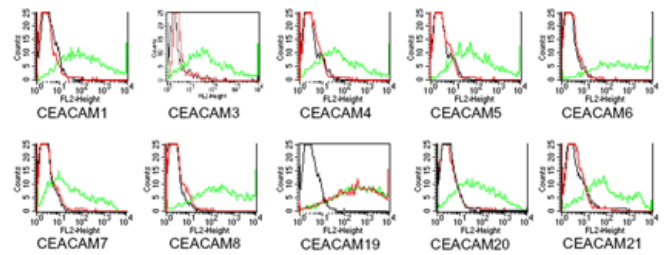
Method: Spectral confocal microscopy of CHO cells using CEACAM19 (human), mAb (HY-8H10) (Prod. No. ALX-805-089). CHO cells were transiently transfected with an expression vector encoding CEACAM19. Binding of CEACAM19 (human), mAb (HY-8H10) was visualized with a FITC-conjugated secondary antibody (green). Actin filaments are labeled with Alexa Fluor-555 Phalloidin (red). Cell nuclei are stained with DAPI (blue).



Specificity testing by flow cytometry. **Method:** FACS analysis of BOSC23 cells using CEACAM19 (human), mAb (HY-8H10) (Prod. No. ALX-805-089). BOSC23 cells were transiently transfected with an expression vector encoding either CEACAM19 (red curve) or an irrelevant protein (control transfectant). Binding of CEACAM19 (human), mAb (HY-8H10) was detected with a PE conjugated secondary antibody. A positive signal was obtained only with CEACAM19 transfected cells.



SDS-PAGE analysis of CEACAM19 (human), mAb (HY-8H10) (Prod. No. ALX-805-089). **Method:** Lane 1: molecular weight marker, Lane 2: 2µg of CEACAM19 (human), mAb (HY-8H10).



Antibody cross-reactivity with members of the CEA family. **Method:** BOSC23 cells were transiently transfected with expression vectors containing either the cDNA of CEACAM1, CEACAM3-8 or CEACAM19-21. Expression of the constructs was tested with monoclonal antibodies known to recognize the corresponding proteins (green curves). An irrelevant monoclonal antibody served as a negative control (black curves). For specificity testing, CEACAM19 (human), mAb (HY-8H10) (Prod. No. ALX-805-089) was tested on all CEACAM transfectants. A positive signal was obtained only with CEACAM19 transfectants (red curve).

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	Carcinoembryonic antigen-related cell adhesion molecule 19, CEAL1
Application	ELISA, Flow Cytometry
Clone	HY-8H10
Crossreactivity	Does not cross-react with human CEACAM1, 3, 4, 5, 6, 7, 8, 20, or 21.
Formulation	Liquid. In PBS, pH 7.2. Contains no preservatives.
Host	Mouse
Immunogen	Vector containing the cDNA of human CEACAM19.
Isotype	IgG1
Purity Detail	Protein G-affinity purified.
Quality Control	Specificity routinely tested by flow cytometry and spectral confocal microscopy on CEACAM19 transfected BOSC23 and CHO cells.
Recommendation Dilutions/Conditions	Flow Cytometry (1.2µg/10 ⁶ cells)Competitive ELISA (1:200-1:400)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Species Reactivity

Human

UniProt ID

Q7Z692

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

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com