

PRODUCT INFORMATION



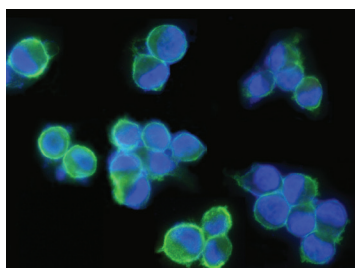
EpCAM Rabbit Monoclonal Antibody (Clone 28)

Item No. 38079

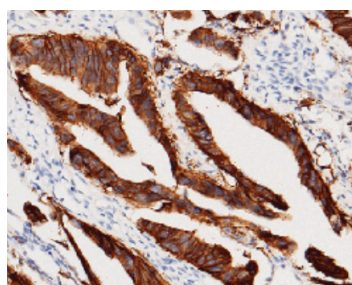
Overview and Properties

Contents:	This vial contains 50, 100 µl, or 1 ml of protein A-affinity purified monoclonal antibody.
Synonyms:	Adenocarcinoma-associated Antigen, CD326, EGP314, Epithelial Cell Adhesion Molecule, Epithelial Cell Surface Antigen, Epithelial Glycoprotein 314, TACSTD1, TROP1, Tumor-associated Calcium Signal Transducer 1
Immunogen:	Recombinant human EpCAM
Cross Reactivity:	(+) EpCAM; (-) CD146, CD171, OBCAM, TROP2
Species Reactivity:	(+) Human
Molecular Weight:	34 kDa
Form:	Liquid
Storage:	-80°C (as supplied)
Stability:	≥1 year
Storage Buffer:	0.2 µm filtered solution in PBS
Clone:	28
Host:	Rabbit
Isotype:	IgG
Applications:	ELISA, ELISA (detection), Flow cytometry (FC), Immunocytochemistry (ICC), Immunofluorescence (IF), Immunohistochemistry paraffin (IHC-P), Immunoprecipitation (IP), and Western blot (WB); the recommended starting dilution is 1:25,000-1:50,000 for ELISA, 1:5,000-1:50,000 for ELISA (detection), 1:100-1:500 for FC, 1:250-1:5,000 for ICC and IF, 1:1,000-1:5,000 for IHC-P, and 1:500-1:1,000 for WB. The recommended starting concentration is 0.2-1 µl/mg for IP. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

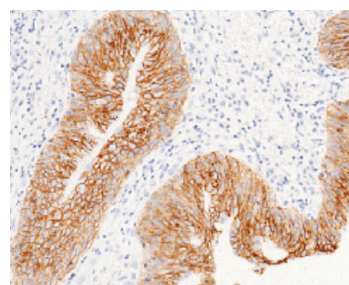
Images



Immunofluorescent staining of human EpCAM in SKBR3 cells. Cells were fixed with 4% paraformaldehyde, blocked with 10% serum, and incubated with EpCAM Rabbit Monoclonal Antibody (Clone 28) at a dilution of 1:500 at 4°C overnight. Cells were stained with the Alexa Fluor® 488-conjugated Goat Anti-Rabbit IgG Secondary Antibody (green) and counterstained with DAPI (blue). Positive staining was localized to the plasma membrane.



Immunohistochemical staining of human EpCAM in formalin-fixed and paraffin-embedded human colon carcinoma using EpCAM Rabbit Monoclonal Antibody (Clone 28) at a dilution of 1:2,500. Positive staining was localized to membrane of colonic gland epithelium.



Immunohistochemical staining of human EpCAM in formalin-fixed and paraffin-embedded human bladder carcinoma using EpCAM Rabbit Monoclonal Antibody (Clone 28) at a dilution of 1:2,500. Positive staining was localized to membrane of transitional epithelium.

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

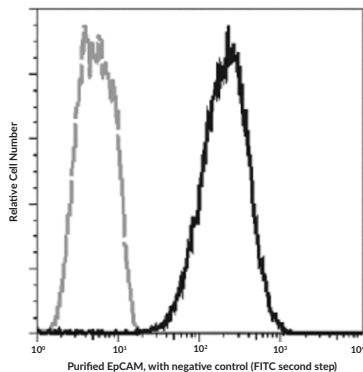
WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/10/2023

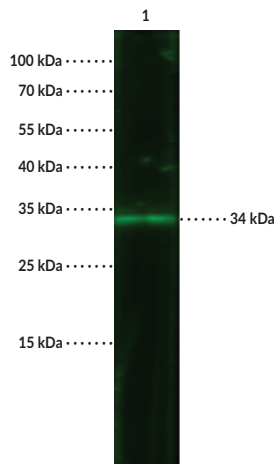
CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION

Images continued

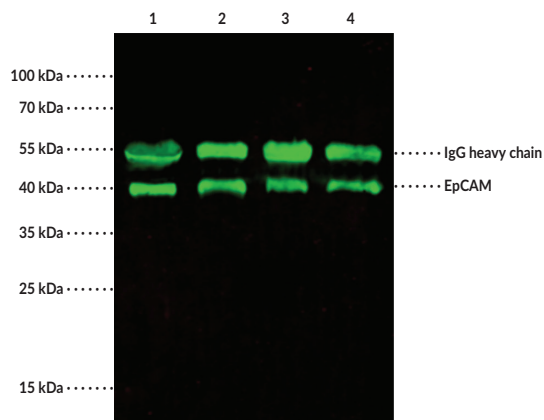


Flow cytometric analysis of EpCAM reactivity on SKBR3 cells. SKBR3 cells were stained with purified EpCAM Rabbit Monoclonal Antibody (Clone 28), followed by a FITC-conjugated second step antibody. The histograms were derived from gated events based on light scattering characteristics of viable cells.



Lane 1: MCF-7 lysate (30 µg)

WB using EpCAM Rabbit Monoclonal Antibody (Clone 28) at a 1:500 dilution.



Lane 1: MCF-7 IP eluate
Lane 2: A431 IP eluate
Lane 3: HepG2 IP eluate
Lane 4: Caco-2 IP eluate

WB using 0.5 µl of EpCAM Rabbit Monoclonal Antibody (Clone 28) and 15 µl of protein G agarose for each 500 µg of four cell lysates.

PRODUCT INFORMATION



Description

Epithelial cell adhesion molecule (EpCAM) is a transmembrane glycoprotein.¹ It is composed of an extracellular domain (EpEX) that contains an N-domain, a thyroglobulin type-IA domain, and a C-domain, a transmembrane domain, and an intracellular domain (EpIC). EpEX undergoes proteolytic cleavage by disintegrin and metalloproteinase domain-containing protein 17 (ADAM17/TACE) to release soluble EpEX, and γ -secretase cleaves EpCAM at two sites to release a soluble extracellular amyloid- β -like fragment and EpIC. EpCAM is ubiquitously expressed in epithelial tissues but is also found in stem and progenitor cells and tight junctions.^{2,3} It is involved in cell adhesion, transport, polarity, proliferation, and motility and signals through novel PKC (nPKC), Wnt, and embryonic RAS (ERAS)/Akt pathways.² Knockdown of *EPCAM* decreases proliferation, migration, and invasion of breast cancer cells *in vitro*.⁴ *Epcam*^{-/-} mice develop congenital tufting enteropathy (CTE), a disease characterized by hemorrhagic diarrhea, colon crypt hyperplasia, intestinal tufts, and villous atrophy.⁵ Increased protein levels of EpCAM are frequently found in tumors from patients with diverse cancer types.⁶ Cayman's EpCAM Rabbit Monoclonal Antibody (Clone 28) can be used for ELISA, ELISA (detection), flow cytometry (FC), immunocytochemistry (ICC), immunofluorescence (IF), immunohistochemistry paraffin (IHC-P), immunoprecipitation (IP), and Western blot (WB) applications. The antibody recognizes EpCAM at 34 kDa from human samples. However, due to glycosylation the antibody may recognize EpCAM at approximately 38-40 kDa.

References

1. Keller, L., Werner, S., and Pantel, K. Biology and clinical relevance of EpCAM. *Cell Stress* **3(6)**, 165-180 (2019).
2. Huang, L., Yang, Y., Yang, F., *et al.* Functions of EpCAM in physiological processes and diseases (Review). *Int. J. Mol. Med.* **42(4)**, 1771-1785 (2018).
3. Ladwein, M., Pape, U.F., Schmidt, D.-S., *et al.* The cell-cell adhesion molecule EpCAM interacts directly with the tight junction protein claudin-7. *Exp. Cell Res.* **309(2)**, 345-357 (2005).
4. Osta, W.A., Chen, Y., Mikhitarian, K., *et al.* EpCAM is overexpressed in breast cancer and is a potential target for breast cancer gene therapy. *Cancer Res.* **64(16)**, 5818-5824 (2004).
5. Guerra, E., Lattanzio, R., La Sorda, R., *et al.* *mTrop1/Epcam* knockout mice develop congenital tufting enteropathy through dysregulation of intestinal E-cadherin/ β -catenin. *PLoS One* **7(11)**, e49302 (2012).
6. Went, P.T., Lugli, A., Meier, S., *et al.* Frequent EpCam protein expression in human carcinomas. *Hum. Pathol.* **35(1)**, 122-128 (2004).