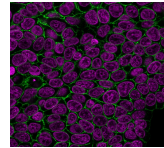


Ep-CAM / CD326 Monoclonal Mouse Antibody (HEA125)



Product Description

Note: All CF® Dye and biotin conjugates of this antibody have been discontinued. The antibody is still available unlabeled, with or without BSA. This antibody recognizes a 40-43 kDa transmembrane epithelial glycoprotein, identified as epithelial specific antigen (ESA), or epithelial cellular adhesion molecule (Ep-CAM). It is expressed on baso-lateral cell surface in most simple epithelia and a vast majority of carcinomas with the exception of adult squamous epithelium, hepatocytes and gastric epithelial cells. Antibody to Ep-CAM has been used to distinguish adenocarcinoma from pleural mesothelioma and hepatocellular carcinoma. It is also useful in distinguishing serous carcinomas of the ovary from mesothelioma. It has been reported that this epithelial antigen plays an important role as a tumor-cell marker in lymph nodes from patients with esophageal carcinoma otherwise classified as node-negative. Primary antibodies are available purified, or with a selection of fluorescent CF® dyes and other labels. CF® dyes offer exceptional brightness and photostability. See the [CF® Dye Brochure](#) for more information. Note: Conjugates of blue fluorescent dyes like CF®405S and CF®405M are not recommended for detecting low abundance targets, because blue dyes have lower fluorescence and can give higher non-specific background than other dye colors. **Stock status:** Because Biotium offers a large number of antibody and conjugation options, primary antibody conjugates may be made to order. Typical lead times are up to one week for CF® dye and biotin conjugates, and up to 2-3 weeks for fluorescent protein and enzyme conjugates. Please email order@biotium.com to inquire about stock status and lead times before placing your order.

References

Note: References for this clone sold by other suppliers may be listed for expected applications.

Br J Cancer (1987) 56: 714-721. (immunoprecipitation; RIA; IHC, frozen)

Product attributes

Antibody number	#0825
Antibody reactivity (target)	CD326, Ep-CAM
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	HEA125
Isotype	IgG2b, kappa
Molecular weight	40-43 kDa
Synonyms	Adenocarcinoma-associated Antigen, Cell Surface Glycoprotein Trop-1, Chromosome 17, ECS-1, EGP2, EGP314, EGP40, Epidermal Surface Molecule, Epithelial Glycoprotein 314, ESA, ESA1, FLOT2, Flotillin-2, KSA, Membrane Component, REG-1, Reggie-1, Reggie-2, Surface Marker-1 (M17S1), TACD1, TROP1, Tumor-associated Calcium Signal Transducer 1 (TACSTD1)
Human gene symbol	TACSTD1
Entrez gene ID	4072
SwissProt	P16422
Unigene	542050
Immunogen	Recombinant human EpCAM protein fragment from the extracellular domain (aa77-202) (exact sequence is proprietary)
Antibody target cellular localization	Exosomes/EVs, Plasma membrane
Species reactivity	Human
Verified antibody applications	Flow (surface) (verified), IF (verified), IHC (FFPE) (verified)
Expected antibody applications	RIA (published for clone), IHC (frozen) (published for clone), IP (published for clone)
Antibody application notes	Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody. Immunohistology (formalin): 0.5-1 ug/mL. Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM citrate buffer, pH 6.0, for 10-20 min followed by cooling at RT for 20 min
Positive control	HT29 cells or Ovarian carcinoma
Shipping condition	Room temperature
Storage Conditions	Store at 2 to 8 °C, Protect fluorescent conjugates from light. Note: store BSA-free antibodies at -10 to -35 °C
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Regulatory status	For research use only (RUO)
Antibody/conjugate formulation	Conjugates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide, HRP conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purified, BSA-free: 1 mg/mL in PBS without azide
Antibody research areas	Cancer, Cell adhesion, Exosomes/EVs
Cell/tissue expression	Epithelial cells
Tumor expression	Adenocarcinoma