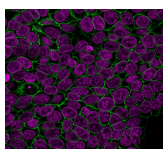


## Ep-CAM / CD326 Monoclonal Mouse Antibody (HEA125)



### 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](mailto:order@biotium.com) to inquire about stock status and lead times before placing your order.

**Catalog number key for antibody number 0825, Anti-CD326|Ep-CAM (HEA125)**

### Product attributes

<b>Antibody number</b>	#0825
<b>Antibody reactivity (target)</b>	CD326, Ep-CAM
<b>Antibody type</b>	Primary
<b>Host species</b>	Mouse
<b>Clonality</b>	Monoclonal
<b>Clone</b>	HEA125
<b>Isotype</b>	IgG2b, kappa
<b>Molecular weight</b>	40-43 kDa
<b>Synonyms</b>	Adenocarcinoma-associated Antigen, Cell Surface Glycoprotein Trop-1, Chromosome 17, ECS-1, EGP2, EGP314, EGP40, Epidermal Surface Antigen 1, Epithelial Cell Adhesion 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)
<b>Human gene symbol</b>	TACSTD1
<b>Entrez gene ID</b>	4072
<b>SwissProt</b>	P16422
<b>Unigene</b>	542050
<b>Immunogen</b>	Recombinant human EpCAM protein fragment from the extracellular domain (aa77-202) (exact sequence is proprietary)
<b>Antibody target cellular localization</b>	Exosomes/EVs, Plasma membrane
<b>Species reactivity</b>	Human
<b>Verified antibody applications</b>	Flow (surface) (verified), IF (verified), IHC (FFPE) (verified)
<b>Expected antibody applications</b>	RIA (published for clone), IHC (frozen) (published for clone), IP (published for clone)
<b>Antibody application notes</b>	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
<b>Positive control</b>	HT29 cells or Ovarian carcinoma
<b>Shipping condition</b>	Room temperature
<b>Storage Conditions</b>	Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C
<b>Shelf life</b>	Guaranteed for at least 24 months from date of receipt when stored as recommended
<b>Regulatory status</b>	For research use only (RUO)
<b>Antibody/conjugate formulation</b>	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
<b>Antibody research areas</b>	Cancer, Cell adhesion, Exosomes/EVs
<b>Cell/tissue expression</b>	Epithelial cells
<b>Product origin</b>	Product may contain either bovine serum albumin (BSA) from bovine serum ( <i>Bos taurus</i> ), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.
<b>Tumor expression</b>	Adenocarcinoma

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel	Dye Features
BNC04	CF@405S	404/431	405	DAPI (microscopy), AF405	<a href="#">CF@405S Features</a>
BNC88	CF@488A	490/515	488	GFP, FITC	<a href="#">CF@488A Features</a>
BNC68	CF@568	562/583	532, 561	RFP, TRITC	<a href="#">CF@568 Features</a>
BNC94	CF@594	593/614	561	Texas Red®	<a href="#">CF@594 Features</a>
BNC40	CF@640R	642/662	633-640	Cy@5	<a href="#">CF@640R Features</a>
BNC47	CF@647	650/665	633-640	Cy@5	<a href="#">CF@647 Features</a>
BNC74	CF@740	742/767	633-685	775/50	<a href="#">CF@740 Features</a>
BNCB	Biotin	N/A	N/A	N/A	
BNUB	Purified	N/A	N/A	N/A	
BNUM	Purified, BSA-free	N/A	N/A	N/A	

Alexa Fluor, Pacific Blue, Pacific Orange, and Texas Red are trademarks or registered trademarks of Thermo Fisher Scientific; Cy is a registered trademark of Cytiva; IRDye, LI-COR, and Odyssey are registered trademarks of LI-COR Bioscience.

## 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)

This datasheet was generated on March 31, 2025 at 12:31:14 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/mouse-monoclonal-anti-ep-cam-cd326-hea125/>