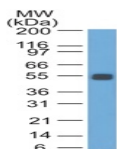




Glowing products for science

# Ep-CAM / CD326 Monoclonal Mouse Antibody (GZ-1)

This item is discontinued.



## Product Description

This item is discontinued. **Catalog number key for antibody number 0381, Anti-CD326|Ep-CAM (GZ-1)**

Call us : [800-304-5357](tel:800-304-5357) Email: [techsupport@biotium.com](mailto:techsupport@biotium.com)

## Product attributes

Antibody number	#0381
Antibody reactivity (target)	CD326, Ep-CAM
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	GZ-1
Isotype	IgG1, kappa
Molecular weight	40-43 kDa
Synonyms	Adenocarcinoma-associated Antigen; Cell Surface Glycoprotein Trop-1; EGP2; EGP314; EGP40; Epithelial Cell Adhesion Molecule; Epithelial Glycoprotein 314; ESA; KSA; TACD1; TACSTD1; TROP1; Tumor-associated Calcium Signal Transducer 1; ECS-1; Epidermal Surface Antigen 1; ESA1; FLOT2; Flotillin-2; M17S1 (Membrane Component; Chromosome 17; Surface Marker-1); REG-1; Reggie-1; Reggie-2
Entrez gene ID	171577
SwissProt	O55159
Immunogen	Rat small intestine cells
Antibody target cellular localization	Plasma membrane
Species reactivity	Rat
Antibody application notes	For coating for ELISA, order Ab without BSA, Optimal dilution and staining procedure for a specific application should be determined by user, Recommended starting concentrations for titration are 1-2 ug/mL for most applications, or 1 ug/million cells/100 uL for flow cytometry
Positive control	Rat small intestine or colon. Rat colon lysate.
Shipping condition	Room temperature
Storage Conditions	Note: store BSA-free antibodies at -10 to -35 °C, Store at 2 to 8 °C, Protect fluorescent conjugates from light
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
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Product origin	Product may contain either bovine serum albumin (BSA) from bovine serum (Bos taurus), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.
Cell/tissue expression	Epithelial cells
Tumor expression	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

Tsubura A et. al. J Cut Pathol, 1992, 19:73-9. |Litvinov SV et. al. J Cell Biol, 1994, 125:437-46. |

This datasheet was generated on August 4, 2025 at 04:14:08 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/ep-cam-cd326-monoclonal-mouse-antibody-gz-1/>