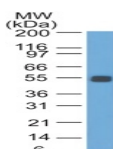


# Ep-CAM / CD326 Monoclonal Mouse Antibody (GZ-1)

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)

## Product attributes

<b>Antibody number</b>	#0381
<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>	GZ-1
<b>Isotype</b>	IgG1, kappa
<b>Molecular weight of antigen</b>	40-43 kDa
<b>Synonyms</b>	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
<b>Entrez gene ID</b>	171577
<b>SwissProt</b>	O55159
<b>Immunogen</b>	Rat small intestine cells
<b>Antibody target cellular localization</b>	Plasma membrane
<b>Species reactivity</b>	Rat
<b>Antibody application notes</b>	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
<b>Positive control</b>	Rat small intestine or colon. Rat colon lysate.
<b>Shipping condition</b>	Room temperature
<b>Storage Conditions</b>	Note: store BSA-free antibodies at -10 to -35 °C, Store at 2 to 8 °C, Protect fluorescent conjugates from light
<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>Shelf life</b>	Guaranteed for at least 24 months from date of receipt when stored as recommended
<b>Product origin</b>	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.
<b>Cell/tissue expression</b>	Epithelial cells
<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

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 June 16, 2026 at 02:00:35 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/>