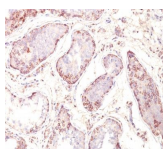


MAGE-A1 Monoclonal Mouse Antibody (MZ2E/838)



Product Description

Recognizes a protein of 42-46 kDa, identified as MAGE-1. This MAb does not cross-react with other members of MAGE-family. Human malignant neoplasms carry rejection antigens that are recognized by the patients' autologous, tumor directed and specific, cytolytic, CD8 T lymphocyte clones (CTL). The MAGE family of genes codes an important group of antigens. It was identified that melanomas and primary glial brain tumors express common melanoma associated antigens (MAAs). Because MAGE-1 is expressed on a significant proportion of human neoplasms of various histological types (melanoma, brain tumors of glial origin, neuroblastoma, non-small cell lung cancer, breast, gastric, colorectal, ovarian, renal cell carcinomas) and not on normal tissues, the encoded antigen may serve as a marker of early detection and target for cancer immunotherapy.

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.

Catalog number key for antibody number 0838, Anti-MAGE-A1 (MZ2E/838)

Product attributes

Antibody number	#0838
Antibody reactivity (target)	MAGE-A1
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	MZ2E/838
Isotype	IgG1, kappa
Molecular weight	42-46 kDa
Synonyms	MZ2 E, cancer/testis antigen 1.1, CT1.1, MAGE1A, MAGEA1, Melanoma antigen family A 1, Melanoma associated antigen 1, Melanoma associated antigen MZ2 E
Human gene symbol	MAGEA1
Entrez gene ID	4100
SwissProt	P43355
Unigene	72879
Immunogen	Recombinant human MAGEA1 protein
Verified antibody applications	IHC (FFPE) (verified)
Antibody target cellular localization	Cytoplasmic
Species reactivity	Human
Antibody application notes	Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Immunofluorescence: 1-2 ug/mL, Immunohistology formalin-fixed 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 minutes, Flow Cytometry 0.5-1 ug/million cells/0.1 mL, Optimal dilution for a specific application should be determined by user
Positive control	Melanoma cell lines. Melanomas, gliomas, neuroblastoma, non-small cell lung cancer, breast, gastric, colorectal, ovarian, and renal cell carcinomas.
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
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.
Tumor expression	Breast cancer, Colorectal cancer, Gastrointestinal cancer, Glioma, Lung cancer, Melanoma, Ovarian cancer

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

This datasheet was generated on November 26, 2024 at 11:42:14 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/monoclonal-anti-mage-a1-mz2e838/>