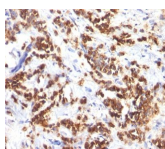


Milk Fat Globulin Monoclonal Mouse Antibody (MFG-06)



Product Description

Recognizes a protein of 40-45 kDa, identified as human milk fat globule membrane protein (HMFG). HMFG is present on normal human breast epithelial cells and cell lines derived from breast carcinomas, as well as to the outer surface of the human milk fat globule. HMFG is considered as a differentiation marker. It is useful as specific breast epithelial marker and can also provide a tool to study the role of the cell surface in normal and neoplastic mammary development.

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 0227, Anti-Milk-Fat-Globulin (MFG-06)

Call us : [800-304-5357](tel:800-304-5357)

Product attributes

Antibody number	#0227
Antibody reactivity (target)	Milk-Fat-Globulin
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	MFG-06
Isotype	IgG1, kappa
Molecular weight of antigen	45 kDa
Synonyms	Breast epithelial antigen BA46; EDIL1; HMFG; Lactadherin; Lactadherin precursor; Medin; MFGE8; MFGM; Milk fat globule-EGF factor 8 (MFG-E8)
Human gene symbol	MFGE8
Entrez gene ID	4240
SwissProt	Q08431
Unigene	3745
Immunogen	Human milk fat globule membrane preparation
Verified antibody applications	IHC (FPPE) (verified)
Antibody target cellular localization	Plasma membrane
Species reactivity	Human
Expected antibody applications	IF (published for clone), WB (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, 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	Normal breast or breast 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
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, Mammary gland

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.

References

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

1. Mol Endocrinol (2010) 24(2): 359-369. (WB)
2. Lab Invest (2014) 94:1260-1272. (IHC, FFPE; IF)
3. Exp Cell Res (2020) 392(1): 112013. (IF)

This datasheet was generated on June 4, 2026 at 11:13:32 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/monoclonal-anti-milk-fat-globulin-mfg-06/>

