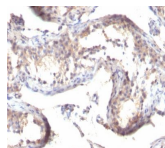


TGF-Alpha Monoclonal Mouse Antibody (MF9)



Product Description

This antibody reacts with the C-terminus of TGF alpha and shows no cross-reaction with EGF and the neuropeptide synenkephalin. The staining with Ab-1 is completely blocked by the peptide used for raising this antibody. TGF-beta (aa50) is a growth factor with 33% homology to EGF, binds to EGFR, activates tyrosine phosphorylation of the receptor, and stimulates cell proliferation. It plays a role in tumor initiation by inducing the reversible transformed phenotype. 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 0644, Anti-TGF-Alpha (MF9)**

Product attributes

Antibody number	#0644
Antibody reactivity (target)	TGF-Alpha
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	MF9
Isotype	IgG1, kappa
Molecular weight	6 kDa
Synonyms	EGF-like TGF; ETGF; TFGA; TGF Type 1; TGFA; Wa1; Waved 1
Human gene symbol	TGFA
Entrez gene ID	7039
SwissProt	P01135
Unigene	170009
Immunogen	A 17-amino acid synthetic peptide from the carboxyl-terminus of rat TGF?
Antibody target cellular localization	Secreted (extracellular)
Verified antibody applications	IHC (FFPE) (verified)
Species reactivity	Cow, Human, Mouse, Rat
Antibody application notes	Immunohistology formalin-fixed 2-4 ug/mL. Staining of formalin-fixed tissues requires boiling tissue sections in 10 mM Tris with 1 mM EDTA, pH 9.0, for 10-20 min followed by cooling at RT for 20 minutes. Flow Cytometry 0.5-1 ug/million cells/0.1 mL. Immunofluorescence 1-2 ug/mL. Optimal dilution for a specific application should be determined by user
Positive control	Jurkat or Ramos cells. Testicular 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
Antibody research areas	Signal transduction

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
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 LI-COR Bioscience.