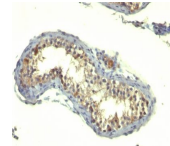


# TGF-Alpha Monoclonal Mouse Antibody (TG86)



## Product Description

This antibody reacts with the TGF alpha and shows no cross-reaction with EGF and the neuropeptide synenkephalin. The staining with this MAb 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](mailto:order@biotium.com) to inquire about stock status and lead times before placing your order.

**Catalog number key for antibody number 0786, Anti-TGF-Alpha (TG86)**

## Product attributes

Antibody number	#0786
Antibody reactivity (target)	TGF-Alpha
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	TG86
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 10-amino acid synthetic peptide (aa34-43; PPVAAAVVSH) from human TGF?
Antibody target cellular	Secreted (extracellular)
Verified antibody applications	IHC (FFPE) (verified)
Species reactivity	Human, Rabbit, Zebrafish
Antibody application notes	Immunohistology formalin-paraffin 2-4 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, 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	<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>
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, of LI-COR Bioscience.