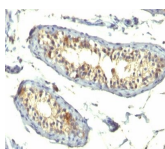


TGF-Alpha Monoclonal Mouse Antibody (TG86 + P/T1)



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.

This antibody is available purified with BSA/azide at 200 ug/mL, or BSA/azide-free at 1 mg/mL.

Catalog number key for antibody number 0790, Anti-TGF-Alpha (TG86 P/T1)

Product attributes

Antibody number	#0790
Antibody reactivity (target)	TGF-Alpha
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	TG86 + P/T1
Isotype	IgM
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? (TG86 & P/T1)
Antibody target cellular localization	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, 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
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.

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 September 15, 2025 at 10:13:19 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/monoclonal-anti-tgfa-alpha-ig86-pt1/>