



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357) Email: [techsupport@biotium.com](mailto:techsupport@biotium.com)

# TGF-Beta Monoclonal Mouse Antibody (TGFB/510)

## Product Description

This MAbs recognizes TGF beta 1, 2 and 3. TGFs inhibit the growth of epithelial cells and stimulate the growth of mesenchymal cells. Most breast lesions, benign and malignant, involve abnormal proliferation and altered architecture of stromal and/or epithelial elements. **Catalog number key for antibody number 0510, Anti-TGF-Beta (TGFB/510)**

## Product attributes

Antibody number	#0510
Antibody reactivity (target)	TGF-Beta
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	TGFB/510
Isotype	IgG1, kappa
Molecular weight	~13 kDa
Synonyms	ARVD; BSC 1 cell growth inhibitor; CED; Cetermin; Diaphyseal dysplasia 1 (DPD1); Glioblastoma derived T cell suppressor factor (G TSF); Latency-associated peptide (LAP); Polyargin; Transforming growth factor beta 1 (TGF beta1 or TGFB1); TGF beta2 (TGFB2); TGF beta3 (TGFB3); TGFbeta (TGFB)
Human gene symbol	TFGB
Entrez gene ID	7040; 7042; 7043
SwissProt	P01137 ; P10600 ; P61812
Unigene	645227
Immunogen	Full length recombinant human TGF-beta2 protein
Antibody target cellular localization	Secreted (extracellular)
Species reactivity	Cow, Dog, Hamster, Human, Monkey, Mouse, Rat
Antibody application notes	For coating for ELISA, order Ab without BSA, Higher concentration may be required for direct detection using primary antibody conjugates than for indirect detection with secondary antibody, Optimal dilution and staining procedure for a specific application should be determined by user, Recommended starting concentrations for titration are 1-2 ug/mL for most applications, or 1 ug/million cells/100 uL for flow cytometry
Positive control	Ovarian carcinoma, lung tumor or skin. PC-3 cell line. Eahy926 cells.
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
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
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
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 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>
BNC74	CF@740	742/767	633-685	775/50	<a href="#">CF@740 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, and Odyssey are registered trademarks of LI-COR Bioscience.

This datasheet was generated on August 28, 2025 at 04:26:32 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/tgf-beta-monoclonal-mouse-antibody-tgfb-510/>