



Glowing products for science

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

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 | 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 December 12, 2025 at 06:55:02 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/tgf-beta-monoclonal-mouse-antibody-tgfb-510/>