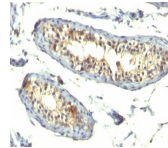


# TGF-Alpha Monoclonal Mouse Antibody (TG86 + P/T1)



## 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.

## Product attributes

<b>Antibody number</b>	#0790
<b>Antibody reactivity (target)</b>	TGF-Alpha
<b>Antibody type</b>	Primary
<b>Host species</b>	Mouse
<b>Clonality</b>	Monoclonal
<b>Clone</b>	TG86 + P/T1
<b>Isotype</b>	IgM
<b>Molecular weight</b>	6 kDa
<b>Synonyms</b>	EGF-like TGF; ETGF; TFGA; TGF Type 1; TGFA; Wa1; Waved 1
<b>Human gene symbol</b>	TGFA
<b>Entrez gene ID</b>	7039
<b>SwissProt</b>	P01135
<b>Unigene</b>	170009
<b>Immunogen</b>	A 10-amino acid synthetic peptide (aa34-43; PPVAAAVVSH) from human TGF $\alpha$ ? (TG86 & P/T1)
<b>Antibody target cellular location</b>	Secreted (extracellular)
<b>Verified antibody applications</b>	IHC (FFPE) (verified)
<b>Species reactivity</b>	Human, Rabbit, Zebrafish
<b>Antibody application notes</b>	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
<b>Positive control</b>	Jurkat or Ramos cells. Testicular Carcinoma
<b>Shipping condition</b>	Room temperature
<b>Storage Conditions</b>	Store at 2 to 8 °C. Note: store BSA-free antibodies at -10 to -35 °C
<b>Shelf life</b>	Guaranteed for at least 24 months from date of receipt when stored as recommended
<b>Regulatory status</b>	For research use only (RUO)
<b>Antibody/conjugate formulation</b>	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
<b>Antibody research areas</b>	Signal transduction