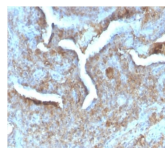


# Thymidine Phosphorylase / PD-ECGF Recombinant Monoclonal Rabbit Antibody (TYMP/2890R)



## Product Description

This antibody recognizes a protein (amino acid 482) of 55 kDa (in vivo 110 kDa homodimer), identified as platelet-derived endothelial growth factor (PD-ECGF), same as thymidine phosphorylase (TP) or gliostatin. In the presence of inorganic orthophosphate, it catalyzes the reversible phospholytic cleavage of thymidine and deoxyuridine to their corresponding bases and 2-deoxyribose-1-phosphate. It is both chemotactic and mitogenic for endothelial cells and a non-heparin binding angiogenic factor present in platelets. Its enzymatic activity is crucial for angiogenic activity (metabolite is angiogenic). Higher levels of serum TP/PD-ECGF are observed in cancer patients. It is also involved in transformation of fluoropyrimidines, cytotoxic agents used in the treatment of a variety of malignancies, into active cytotoxic metabolites (e.g. 5'-deoxy-5-fluorouridine to 5-FU). High intra-cellular levels of TP/PD-ECGF are associated with increased chemosensitivity to such antimetabolites

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.

## Product attributes

Antibody number	#2890
Antibody reactivity (target)	PD-ECGF, Thymidine Phosphorylase
Antibody type	Primary
Host species	Rabbit
Clonality	Recombinant Monoclonal
Clone	TYMP/2890R
Isotype	IgG
Molecular weight	55 kDa
Synonyms	ECGF; ECGF1; Gliostatin; hPD-ECGF; MEDPS1; MNGIE; MTDPS1; PD-ECGF; PDECGF; Platelet-derived endothelial cell growth factor; TdRPase; Thymidine phosphorylase; TP; Tymp
Human gene symbol	TYMP
Entrez gene ID	1890
SwissProt	P19971
Unigene	180903
Immunogen	Recombinant full-length human Thymidine Phosphorylase (TP / PD-ECGF) protein
Verified antibody applications	IHC (FPPE) (verified)
Antibody target cellular localization	Nucleus & cytoplasm
Species reactivity	Human, Mouse, Rat
Positive control	HUVEC cells, Breast, Bladder, Lung or Kaposi tumors
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	Angiogenesis, Metabolism

This datasheet was generated on September 2, 2024 at 04:20:19 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/thymidine-phosphorylase-pd-ecgf-recombinant-monoclonal-rabbit-antibody-tymp-2890r/>