

PRODUCT INFORMATION



TP Receptor (mouse) Polyclonal Antibody

Item No. 101882

Overview and Properties

Contents:	This vial contains 500 µl of peptide affinity-purified polyclonal antibody.
Synonyms:	Prostanoid TP Receptor, Thromboxane A ₂ Receptor, TXA ₂ Receptor
Immunogen:	Synthetic peptide from an internal region of the mouse TP receptor
Cross Reactivity:	(+) TP _α , TP _β
Species Reactivity:	(+) Human, bovine, monkey, mouse, rat
Uniprot No.:	P30987
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥3 years
Storage Buffer:	TBS, pH 7.4, with 50% glycerol, 0.1% BSA, and 0.02% sodium azide
Host:	Rabbit
Application:	Western blot (WB); the recommended starting dilution is 1:200. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Description

The TP receptor is a G protein-coupled receptor that mediates the action of thromboxane A₂ (TXA₂).¹ Alternative splicing of TP pre-mRNA produces two receptor variants, TP_α and TP_β, which differ in sequence and the length of their carboxy terminal tail. TP receptors are expressed primarily in platelets but are also expressed in placenta, vascular smooth muscle, brain, small intestine, thymus, kidney, lung, spleen, and monocytes.¹⁻⁴ It signals in a cell type- and expression-dependent manner and has roles in cardiovascular and hemodynamic function.^{1,5} *Tp*^{-/-} mice exhibit coagulation defects, altered vascular responses to TXA₂ and arachidonic acid, and increased susceptibility to arachidonic acid-induced shock.⁶ Cayman's TP Receptor (mouse) Polyclonal Antibody can be used for Western blot (WB). The antibody recognizes the TP_α and TP_β receptors from human, bovine, monkey, mouse, and rat samples. The apparent molecular weight for TP receptors has been reported from 37 kDa to 70 kDa, depending on different degrees of glycosylation.⁷⁻⁹

References

1. Santilli, F., Mucci, L., and Davi, G. *Intern. Emerg. Med.* **6**(3), 203-212 (2011).
2. Namba, T., Sugimoto, Y., Hirata, M., et al. *Biochem. Biophys. Res. Commun.* **184**(3), 1197-1203 (1992).
3. Båtshake, B., Nilsson, C., and Sundelin, J. *Biochem. Biophys. Res. Commun.* **256**(2), 391-397 (1999).
4. Mais, D.E., True, T.A., and Martinelli, M.J. *Eur. J. Pharmacol.* **227**(3), 267-274 (1992).
5. Petri, M.H., Tellier, C., Michiels, C., et al. *Biochem. Biophys. Res. Commun.* **441**(2), 393-398 (2013).
6. Thomas, D.W., Mannon, R.B., Mannon, P.J., et al. *J. Clin. Invest.* **102**(11), 1994-2001 (1998).
7. Wang, G.-R., Zhu, Y., Halushka, P.V., et al. *Proc. Natl. Acad. Sci. USA* **95**(9), 4888-4893 (1998).
8. Takahashi, N., Takeuchi, K., Abe, T., et al. *Endocrinology* **137**(11), 5170-5173 (1996).
9. Habib, A., Fitzgerald, G.A., and Maclouf, J. *J. Biol. Chem.* **274**(5), 2645-2651 (1999).

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/02/2023

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM