

PRODUCT INFORMATION



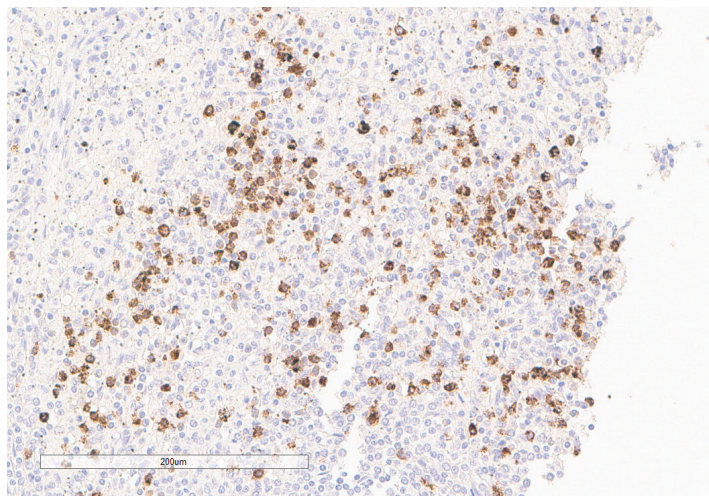
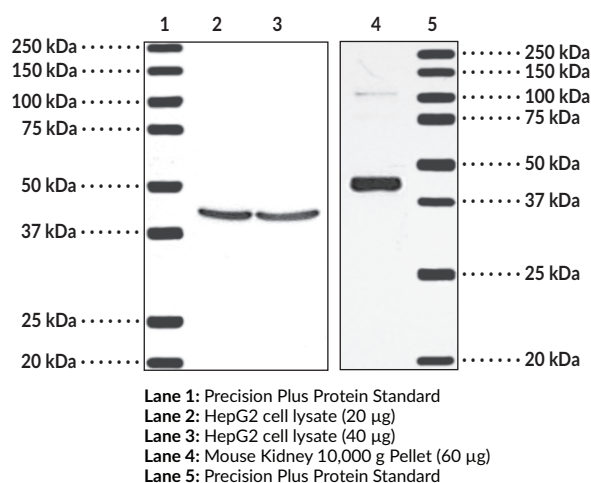
TP Receptor (human) Polyclonal Antibody

Item No. 10004452

Overview and Properties

Contents:	This vial contains peptide affinity-purified IgG.
Synonyms:	Thromboxane A ₂ Receptor, TXA ₂ Receptor
Immunogen:	Synthetic peptide from the C-terminal region of human TP receptor
Species Reactivity:	(+) Human, COS-7 (African green monkey), mouse, and rat TP receptor
Uniprot No.:	P21731
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥3 years
Storage Buffer:	PBS, pH 7.2, with 50% glycerol and 0.02% sodium azide
Host:	Rabbit
Applications:	Immunocytochemistry (ICC), Immunohistochemistry (IHC), and Western blot (WB); the recommended starting dilution is 1:200. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Images



Immunohistochemistry analysis of formalin-fixed, paraffin-embedded (FFPE) human liver tissue after heat induced antigen retrieval in pH 6.0 citrate buffer. After incubation with TP Receptor (human) Polyclonal Antibody (Item No. 10004452), at a 1:200 dilution, slides were incubated with biotinylated secondary antibody, followed by alkaline phosphatase-streptavidin and chromogen (DAB).

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, 10/11/2023

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

PRODUCT INFORMATION



Description

Thromboxane A₂ (TXA₂) is a potent vasoconstrictor and activator of platelet aggregation. TXA₂ elicits its effects via a 7-transmembrane domain G protein-coupled receptor, the TP receptor.¹ The TP receptor is highly expressed in platelets and is relatively less abundant in tissues such as lung, kidney, brain, spleen, thymus, monocytes, uterus, and placenta.²⁻⁵ The apparent molecular weight for TP receptors has been reported from 37 kDa to 70 kDa, depending on different degrees of glycosylation.⁵⁻⁷ However Cayman's antibody consistently detects the TP receptor at 55 and 64 kDa in platelet, kidney, and Cos-7 cell samples. Cayman Chemical's TP receptor polyclonal antibody can be used for western blot and immunocytochemical analysis for both the TPA and TPB splice variants of the receptor on samples of human, murine, rat, and COS-7 (African green monkey) origin.

References

1. Hirata, M., Hayashi, Y., Ushikubi, F., *et al.* Cloning and expression of cDNA for a human thromboxane A₂ receptor. *Nature* **349**(6310), 617-620 (1991).
2. Namba, T., Sugimoto, Y., Hirata, M., *et al.* Mouse thromboxane A₂ receptor: cDNA cloning, expression and Northern blot analysis. *Biochem. Biophys. Res. Commun.* **184**(3), 1197-1203 (1992).
3. Båttshake, B., Nilsson, C., and Sundelin, J. Structure and expression of the murine thromboxane A₂ receptor gene. *Biochem. Biophys. Res. Commun.* **256**(2), 391-397 (1999).
4. Mais, D.E., True, T.A., and Martinelli, M.J. Characterization by photoaffinity labelling of the human platelet thromboxane A₂/prostaglandin H₂ receptor: Evidence for N-linked glycosylation. *Eur. J. Pharmacol.* **227**(3), 267-274 (1992).
5. Takahashi, N., Takeuchi, K., Abe, T., *et al.* Immunolocalization of rat thromboxane receptor in the kidney. *Endocrinology* **137**(11), 5170-5173 (1996).
6. Wang, G.-R., Zhu, Y., Halushka, P.V., *et al.* Mechanism of platelet inhibition by nitric oxide: *In vivo* phosphorylation of thromboxane receptor by cyclic GMP-dependent protein kinase. *Proc. Natl. Acad. Sci. USA* **95**(9), 4888-4893 (1998).
7. Habib, A., Fitzgerald, G.A., and Maclouf, J. Phosphorylation of the thromboxane receptor α , the predominant isoform expressed in human platelets. *J. Biol. Chem.* **274**(5), 2645-2651 (1999).

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