

# PRODUCT INFORMATION



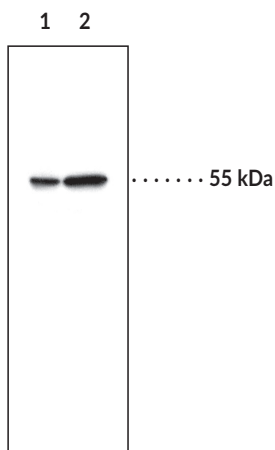
## Thromboxane Synthase Polyclonal Antibody

Item No. 160715

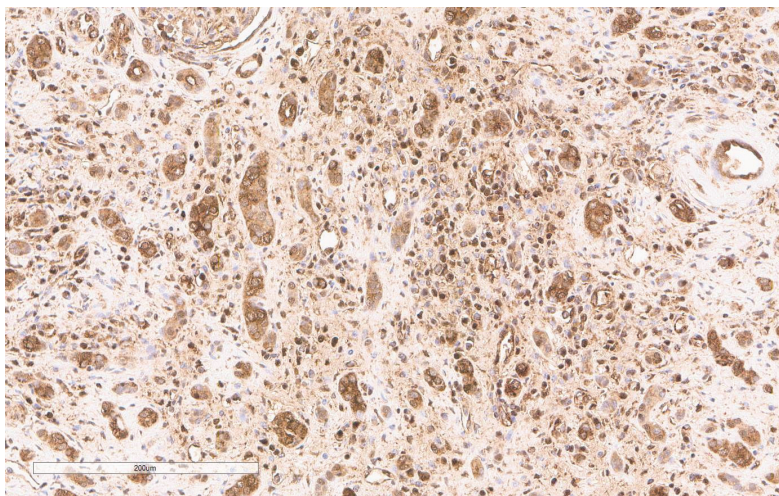
### Overview and Properties

<b>Contents:</b>	This vial contains 500 µl of peptide affinity-purified polyclonal antibody.
<b>Synonyms:</b>	Cytochrome P450 5A1, TBXAS1, Thromboxane A Synthase, TXAS, TXS, TX Synthase, TXA Synthase
<b>Immunogen:</b>	Synthetic peptide from the internal region of human TX synthase
<b>Species Reactivity:</b>	(+) Human; other species not tested
<b>Uniprot No.:</b>	P24557
<b>Form:</b>	Liquid
<b>Storage:</b>	-20°C (as supplied)
<b>Stability:</b>	≥3 years
<b>Storage Buffer:</b>	PBS, pH 7.2, with 50% glycerol and 0.02% sodium azide
<b>Host:</b>	Rabbit
<b>Applications:</b>	Immunohistochemistry (IHC) and Western blot (WB); the recommended starting dilution for IHC is 1:40 and 1:200 for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

### Images



Lane 1: Human platelet microsomes (10 µg)  
Lane 2: Human platelet microsomes (30 µg)



Immunohistochemistry analysis of formalin-fixed, paraffin-embedded (FFPE) human kidney tissue after heat induced antigen retrieval in pH 6.0 citrate buffer. After incubation with Thromboxane Synthase Polyclonal Antibody (Item No. 160715) at a 1:40 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, 11/16/2023

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 (TX) synthase catalyzes the conversion of prostaglandin H<sub>2</sub> to TXA<sub>2</sub>, which is a potent vasoconstrictor and inducer of platelet aggregation. The enzyme has been cloned from a variety of species including human, pig, rat, and mouse and has a molecular weight of approximately 60,000.<sup>1-5</sup> Northern blot analysis reveals that the gene is expressed in lung, kidney and spleen.<sup>4,5</sup>

## References

---

1. Ohashi, K., Ruan, K.-H., Kulmacz, R.J., *et al.* Primary structure of human thromboxane synthase determined from the cDNA sequence. *J. Biol. Chem.* **267**(2), 789-793 (1992).
2. Yokoyama, C., Miyata, A., Ihara, H., *et al.* Molecular cloning of human platelet thromboxane A synthase. *Biochem. Biophys. Res. Commun.* **178**(3), 1479-1484 (1991).
3. Shen, R.-F., Zhang, L., Baek, S.J., *et al.* The porcine thromboxane synthase-encoding cDNA: Sequence, mRNA expression and enzyme production in Sf9 insect cells. *Gene* **140**(2), 261-265 (1994).
4. Tone, Y., Miyata, A., Hara, S., *et al.* Abundant expression of thromboxane synthase in rat macrophages. *FEBS Lett.* **340**(3), 241-244 (1994).
5. Zhang, L., Chase, M.B., and Shen, R.F. Molecular cloning and expression of murine thromboxane synthase. *Biochem. Biophys. Res. Commun.* **194**(2), 741-748 (1997).

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