

PRODUCT INFORMATION



Thromboxane A Synthase (human, recombinant)

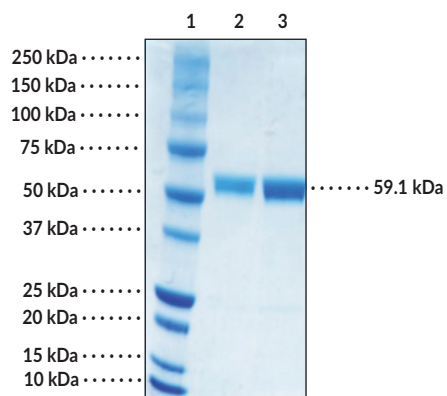
Item No. 33678

Overview and Properties

Synonyms: CYP5A1, Cytochrome P450 5A1, TBXAS1, TXAS, TXA Synthase
Source: Active recombinant C-terminal His-tagged thromboxane A synthase expressed in *E. coli*
Amino Acids: 30-533 with a modified N-terminus and F-G loop
Uniprot No.: P24557
Molecular Weight: 59.1 kDa
Storage: -80°C (as supplied)
Stability: ≥6 months
Purity: *batch specific* (≥90% estimated by SDS-PAGE)
Supplied in: 50 mM Tris, pH 7.4, with 5% glycerol, 0.2 mM DTT, and 0.2 mM EDTA
Protein
Concentration: *batch specific* mg/ml
Activity: Confirmed

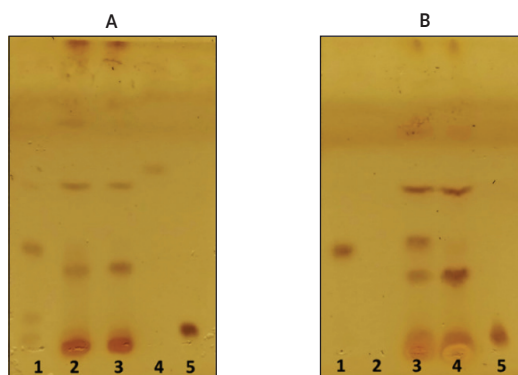
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Images



Lane 1: MW Markers
Lane 2: Thromboxane A Synthase (2 µg)
Lane 3: Thromboxane A Synthase (4 µg)

SDS-PAGE Analysis of Thromboxane A Synthase.



Panel A: Assay using 18 µM TXAS

Panel B: Assay using 9 µM TXAS

Lane 1: PGH₂
Lane 2: PGH₂ conversion for 30 min.
Lane 3: PGH₂ conversion for one hour.
Lane 4: 12(S)-HHTrE (Item No. 34590)
Lane 5: TXB₂ (Item No. 19030)

Lane 1: PGH₂
Lane 2: NO substrate control
Lane 3: PGH₂ conversion for 30 min.
Lane 4: PGH₂ conversion for one hour.
Lane 5: TXB₂ (Item No. 19030)

Thin layer chromatographic analysis of Prostaglandin H₂ (PGH₂; Item No. 17020) conversion by recombinant TXAS (human, recombinant). *In vitro* conversion of PGH₂ (50 µM) was done using 18 µM (Panel A) and 9 µM (Panel B) in 250 µL of assay buffer (10 mM potassium phosphate buffer, pH 7.4, with 0.2 mM DTT, 0.5 mM EDTA and 10% glycerol). The concentration of the protein was based on the Ferrous-CO complexed form of the protein. Analysis of the product was performed in a TLC plate using solvent 40:60:1 ethyl acetate:heptane:acetic acid at 4°C. The plates were visualized by spraying 5% sulfuric acid in methanol.

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, 02/16/2023

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 A synthase (TXAS), also known as cytochrome P450 (CYP) isoform CYP5A1, is the enzyme that catalyzes the isomerization of prostaglandin H₂ (PGH₂; Item No. 17020) into thromboxane A₂ (TXA₂), a vasoconstrictor and an inducer of platelet aggregation.^{1,2} It also catalyzes the cleavage of PGH₂ into malondialdehyde (MDA) and 12(S)-hydroxyheptadecatrienoic acid (12(S)-HHT; Item No. 34590), a leukotriene B₄ (LTB₄) receptor 2 (BLT₂) agonist that also has a role in platelet aggregation.¹ TXAS exists as a monomer and is composed of an N-terminal membrane anchor domain, a heme-binding catalytic residue, and several substrate-binding residues.³⁻⁵ It is expressed in numerous cells, including platelets, monocytes, and macrophages, as well as several tissues, and is localized to the endoplasmic reticulum.³ TXAS undergoes suicide inactivation during catalysis. Mice deficient in TXAS exhibit prolonged bleeding time and defective platelet aggregation.⁶ Cayman's Thromboxane A Synthase (human, recombinant) protein can be used for enzyme activity assay and Western blot (WB) applications. To construct this protein, the N-terminal amino acids 1-29 of human CYP5A1 were removed and replaced with a hydrophilic sequence. The F-G loop of TXAS was modified using sequences from rabbit CYP2C5 and CYP2C3. The C-terminal of this protein contains 6x-His tag followed by a stop codon.

References

1. Okuno, T. and Yokomizo, T. Biological functions of 12(S)-hydroxyheptadecatrienoic acid as a ligand of leukotriene B₄ receptor 2. *Inflamm. Regen.* **38**, 29 (2018).
2. Hajeyah, A.A., Griffiths, W.J., Wang, Y., et al. The biosynthesis of enzymatically oxidized lipids. *Front. Endocrinol. (Lausanne)* **11**, 591819 (2020).
3. Ershov, P.V., Yablokov, E., Zgoda, V., et al. A new insight into subinteractomes of functional antagonists: Thromboxane (CYP5A1) and prostacyclin (CYP8A1) synthases. *Cell Biol. Int.* **45(6)**, 1175-1182 (2021).
4. Ruan, K.H., Milfield, K., Kulmacz, R.J., et al. Comparison of the construction of a 3-D model for human thromboxane synthase using P450cam and BM-3 as templates: Implications for the substrate binding pocket. *Protein Eng.* **7(11)**, 1345-1351 (1994).
5. Seo, M.-J. and Oh, D.-K. Prostaglandin synthases: Molecular characterization and involvement in prostaglandin biosynthesis. *Prog. Lipid Res.* **66**, 50-68 (2017).
6. Yu, I.-S., Lin, S.-R., Huang, C.-C., et al. TXAS-deleted mice exhibit normal thrombopoiesis, defective hemostasis, and resistance to arachidonate-induced death. *Blood* **104(1)**, 135-142 (2004).

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