

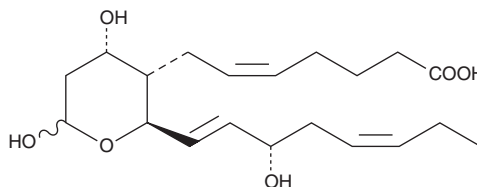
PRODUCT INFORMATION



Thromboxane B₃

Item No. 19990

CAS Registry No.: 71953-80-5
Formal Name: 9 α ,11,15S-trihydroxy-thromba-5Z,13E,17Z-trien-1-oic acid
Synonyms: Δ^{17} -TXB₂, TXB₃
MF: C₂₀H₃₂O₆
FW: 368.5
Purity: $\geq 95\%$
Supplied as: A solution in methyl acetate
Storage: -20°C
Stability: ≥ 2 years



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

Laboratory Procedures

Thromboxane B₃ (TXB₃) is supplied as a solution in methyl acetate. To change the solvent, simply evaporate the methyl acetate under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of TXB₃ in these solvents is approximately 100, 25, and 50 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. If an organic solvent-free solution of TXB₃ is needed, it can be prepared by evaporating the methyl acetate and directly dissolving the neat oil in aqueous buffers. The solubility of TXB₃ in PBS, pH 7.2, is approximately 100 μ g/ml. We do not recommend storing the aqueous solution for more than one day.

Description

TXB₃ is the stable hydrolysis product of TXA₃ synthesized from eicosapentaenoic acid (EPA; Item No. 90110) by COX and thromboxane synthase. It is biosynthesized in various tissues such as seminal vesicles, lung, PMNL, and ocular tissues.^{1,2}

References

1. Kulkarni, P.S., Kaufman, P.L., Srinivasan, B.D. Eicosapentaenoic acid metabolism in cynomolgus and rhesus conjunctiva and eyelid. *J. Ocul. Pharmacol.* **3**, 349-356 (1987).
2. Kulkarni, P.S. and Srinivasan, B.D. Eicosapentaenoic acid metabolism in human and rabbit anterior uvea. *Prostaglandins* **31**, 1159-1164 (1986).

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, 06/30/2017

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