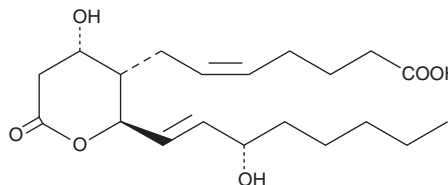


PRODUCT INFORMATION



11-dehydro Thromboxane B₂ Item No. 19500

CAS Registry No.: 67910-12-7
Synonyms: 11-dehydro TXB₂, 11-keto TXB₂
Formal Name: 9 α ,15S-dihydroxy-11-oxothromba-5Z,13E-dien-1-oic acid
MF: C₂₀H₃₂O₆
FW: 368.5
Purity: \geq 98%
Supplied as: A solution in methyl acetate
Storage: -20°C
Stability: \geq 2 years



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

Laboratory Procedures

11-dehydro TXB₂ 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 11-dehydro TXB₂ in these solvents is approximately 100, 25, 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 11-dehydro TXB₂ is needed, it can be prepared by evaporating the methyl acetate and directly dissolving the neat oil in aqueous buffers. The solubility of 11-dehydro TXB₂ in PBS (pH 7.2) is approximately 100 μ g/ml. Avoid adding 11-dehydro TXB₂ to basic solutions (pH 7.4), since base treatment will cause the hydrolysis of the lactone function of 11-dehydro TXB₂. We do not recommend storing the aqueous solution for more than one day.

Description

TXB₂ is released in substantial quantities from aggregating platelets and metabolized during circulation to 11-dehydro TXB₂ and 2,3-dinor TXB₂.¹ 11-dehydro TXB₂ is one of the main plasma metabolites of TXB₂ and can be used as a marker for *in vivo* TXA₂ synthesis.¹⁻⁴ The mean plasma level of 11-dehydro TXB₂ in human males is 0.9-4.3 pg/ml and the half-life is 45-60 minutes.²⁻⁴ Urinary concentrations of 11-dehydro TXB₂ are approximately 30-70 ng/mmol creatinine.^{5,6}

References

1. Ciabattini, G., Pugliese, F., Davi, G., *et al. Biochim. Biophys. Acta* **992**, 66-70 (1989).
2. Fitzgerald, G.A., Lawson, J., Blair, I.A., *et al. Adv. Prostaglandin Thromboxane Leukotriene Res.* **15**, 87-90 (1985).
3. Takasaki, W., Nakagawa, A., Tanaka, Y., *et al. Thromb. Res.* **63**, 331-341 (1991).
4. Catella, F., Healy, D., Lawson, J.A., *et al. Proc. Natl. Acad. Sci. USA* **83**, 5861-5865 (1986).
5. Lellouche, F., Fradin, A., FitzGerald, G., *et al. Prostaglandins* **40**, 297-310 (1990).
6. Perneby, C., Granstrom, E., Beck, O., *et al. Thromb. Res.* **96**, 427-436 (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, 11/13/2023

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