

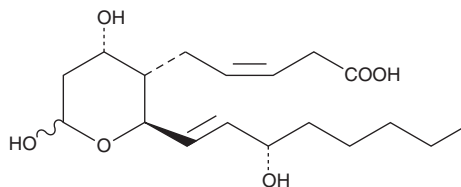
PRODUCT INFORMATION



2,3-dinor Thromboxane B₂

Item No. 19050

CAS Registry No.: 63250-09-9
Formal Name: 9 α ,11,15S-trihydroxy-2,3-dinor-thromba-5Z,13E-dien-1-oic acid
Synonym: 2,3-dinor TXB₂
MF: C₁₈H₃₀O₆
FW: 342.4
Purity: $\geq 95\%$
Supplied as: A 100 $\mu\text{g/ml}$ solution in methyl acetate
Storage: -20°C
Stability: ≥ 4 years



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

Laboratory Procedures

2,3-dinor Thromboxane B₂ (2,3-dinor 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 2,3-dinor TXB₂ 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 2,3-dinor TXB₂ is needed, it can be prepared by evaporating the methyl acetate and directly dissolving the neat oil in aqueous buffers. The solubility of 2,3-dinor TXB₂ in PBS (pH 7.2) is approximately 100 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Thromboxane B₂ (TXB₂) is released in substantial quantities from aggregating platelets and metabolized during circulation to 11-dehydro TXB₂ and 2,3-dinor TXB₂.¹ 2,3-dinor TXB₂ is an abundant urinary metabolite of TXB₂ and can be used as a marker for *in vivo* TXA₂ synthesis.¹⁻³ In healthy male volunteers, the median excretion of 2,3-dinor TXB₂ is 10.3 ng/hour (138 pg/mg creatine).²

References

1. Ciabattoni, G., Pugliese, F., Davi, G., *et al.* Fractional conversion of thromboxane B₂ to urinary 11-dehydrothromboxane B₂ in man. *Biochim. Biophys. Acta* **992**(1), 66-70 (1989).
2. Lawson, J.A., Brash, A.R., Doran, J., *et al.* Measurement of urinary 2,3-dinor-thromboxane B₂ and thromboxane B₂ using bonded-phase phenylboronic acid columns and capillary gas chromatography-negative-ion chemical ionization mass spectrometry. *Anal. Biochem.* **150**(2), 463-470 (1985).
3. Lawson, J.A., Patrono, C., Ciabattoni, G., *et al.* Long-lived enzymatic metabolites of thromboxane B₂ in the human circulation. *Anal. Biochem.* **155**(1), 198-205 (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, 05/14/2026

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