

PRODUCT INFORMATION

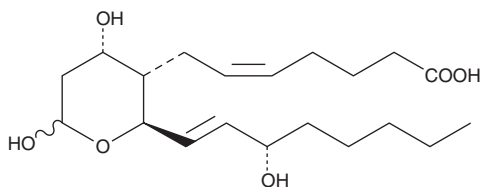


Thromboxane B₂ Quant-PAK

Item No. 10006832

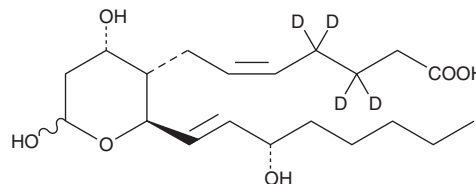
Thromboxane B₂

CAS Registry No.: 54397-85-2
Formal Name: 9 α ,11,15S-trihydroxythromba-5Z,13E-dien-1-oic acid
MF: C₂₀H₃₄O₆
FW: 370.5
Purity: \geq 98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 2 years



Thromboxane B₂-d₄

CAS Registry No.: 1346112-79-5
Formal Name: (5Z)-7-[[[(2R,3S,4S,6R)-tetrahydro-4,6-dihydroxy-2-[(1E,3S)-3-hydroxy-1-octen-1-yl]-2H-pyran-3-yl]-5-heptenoic-3,3,4,4-d₄ acid
MF: C₂₀H₃₀D₄O₆
FW: 374.5
Chemical Purity: \geq 98%
Deuterium Incorporation: \geq 99% deuterated forms (d₁-d₄); \leq 1% d₀
Supplied as: A solution in methyl acetate
Storage: -20°C
Stability: \geq 1 year



Laboratory Procedures

This thromboxane B₂ (TXB₂) Quant-PAK contains 50 μ g of TXB₂-d₄ and 2-4 mg of TXB₂ (please see the vial for exact amount and concentration).

TXB₂ is supplied as a crystalline solid. A stock solution may be made by dissolving the TXB₂ in an organic solvent purged with an inert gas. TXB₂ is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of TXB₂ in these solvents is approximately 100, 25, and 50 mg/ml, respectively.

TXB₂-d₄ 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 TXB₂-d₄ in these solvents is approximately 100, 25, and 50 mg/ml, respectively.

TXB₂-d₄ contains four deuterium atoms at the 3, 3, 4, and 4 positions. It is intended for use as an internal standard for the quantification of TXB₂ by GC- or LC-MS. The accuracy of the sample weight in the TXB₂-d₄ vial is between 5% over and 2% under the weight indicated on the vial. For better precision we have provided a precisely weighed unlabeled TXB₂, with the precise weight (2-4 mg) indicated on the vial. Using this vial the deuterated standard can be quantified by constructing a standard curve of peak intensity ratios (deuterated *versus* unlabeled).

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, 11/30/2021

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