

PRODUCT INFORMATION



Thromboxane B₂-biotin

Item No. 9002298

Formal Name: (Z)-7-((2R,3S,4S)-4,6-dihydroxy-2-((S,E)-3-hydroxyoct-1-en-1-yl)tetrahydro-2H-pyran-3-yl)-N-(5-(5-((3aS,4S,6aR)-2-oxohexahydro-1H-thieno[3,4-d]imidazol-4-yl)pentanamido)pentyl)hept-5-enamide

Synonym: TXB₂-biotin

MF: C₃₅H₆₀N₄O₇S

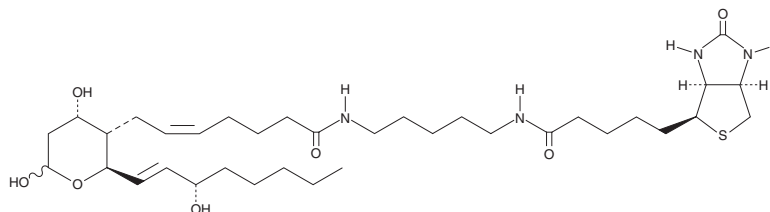
FW: 681.0

Purity: ≥90%

Supplied as: A solution in ethanol

Storage: -20°C

Stability: ≥2 years



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

Description

Thromboxane A₂ (TXA₂) is produced from arachidonic acid (Item No. 90010) by many cells and causes irreversible platelet aggregation and contraction of vascular and bronchial smooth muscle.¹⁻³ TXA₂ is rapidly hydrolyzed non-enzymatically to TXB₂ (Item No. 19030), which is then quickly metabolized to urinary metabolites for clearance by the kidneys. TXB₂-biotin is an affinity probe which allows TXB₂ to be detected through an interaction with the biotin ligand. It was designed to allow TXB₂ to be detected in complexes with protein binding partners.

References

1. Hamberg, M., Svensson, J., and Samuelsson, B. Thromboxanes: A new group of biologically active compounds derived from prostaglandin endoperoxides. *Proc. Natl. Acad. Sci. USA* **72(8)**, 2994-2998 (1975).
2. Ellis, E.F., Oelz, O., Roberts, L.J., II, *et al.* Coronary arterial smooth muscle contraction by a substance released from platelets: Evidence that it is thromboxane A₂. *Science* **193(4258)**, 1135-1137 (1976).
3. Salzman, P.M., Salmon, J.A., and Moncada, S. Prostacyclin and thromboxane A₂ synthesis by rabbit pulmonary artery. *J. Pharmacol. Exp. Ther.* **215(1)**, 240-247 (1980).

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, 02/21/2024

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