

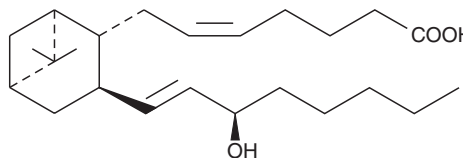
# PRODUCT INFORMATION



## 15(R)-Pinane Thromboxane A<sub>2</sub>

Item No. 10008510

**CAS Registry No.:** 71154-83-1  
**Formal Name:** (5Z)-7-[(1S,2R,3R,5S)-3-[(1E,3R)-3-hydroxy-1-octen-1-yl]-6,6-dimethylbicyclo[3.1.1]hept-2-yl]-5-heptenoic acid  
**Synonyms:** 15-*epi*-Pinane Thromboxane A<sub>2</sub>, 15-*iso*-Pinane Thromboxane A<sub>2</sub>, 15(R)-PTA<sub>2</sub>, 15-*epi*-PTA<sub>2</sub>, 15-*iso*-PTA<sub>2</sub>  
**MF:** C<sub>24</sub>H<sub>40</sub>O<sub>3</sub>  
**FW:** 376.6  
**Purity:** ≥98%  
**Supplied as:** A solution in ethanol  
**Storage:** -20°C  
**Stability:** ≥1 year



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

### Laboratory Procedures

15(R)-Pinane thromboxane A<sub>2</sub> (15(R)-PTA<sub>2</sub>) is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. Solvents such as DMSO and dimethyl formamide purged with an inert gas can be used. The solubility of 15(R)-PTA<sub>2</sub> in these solvents is approximately 25 and 50 mg/ml, respectively.

If an organic solvent-free solution of 15(R)-PTA<sub>2</sub> is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. The solubility of 15(R)-PTA<sub>2</sub> in PBS, pH 7.2, is approximately 0.15 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

15(R)-PTA<sub>2</sub> is the (R)-epimer of pinane thromboxane A<sub>2</sub> (PTA<sub>2</sub>; Item No. 19020). 15(R)-PTA<sub>2</sub> does not inhibit collagen-induced platelet aggregation (IC<sub>50</sub> = 120-130 μM).<sup>1</sup> It does not affect gastric tone in isolated rat gastric fundus when used at concentrations of 0.5 or 1.5 μg/ml and is less effective than PTA<sub>2</sub> at inhibiting prostaglandin-induced contraction of isolated rat stomach muscle.<sup>2</sup>

### References

1. Caton, M.P.L., Stuttle, K.A.J., Tuffin, D.P., *et al.* Comparative action of carbocyclic thromboxane A<sub>2</sub> stereoisomers on platelets. *Eur. J. Med. Chem.* **35(12)**, 1099-1107 (2000).
2. Bennett, A. and Sanger, G.J. Pinane thromboxane A<sub>2</sub> analogues are non-selective prostanoid antagonists in rat and human stomach muscle. *Br. J. Pharmacol.* **77(4)**, 591-596 (1982).

#### 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, 1/03/2018

#### 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