

PRODUCT INFORMATION



I-BOP

Item No. 19600

CAS Registry No: 128719-90-4
Formal Name: (5Z)-7-[(1S,2R,3R,4R)-3-[(1E,3R)-3-hydroxy-4-(4-iodophenoxy)-1-buten-1-yl]-7-oxabicyclo[2.2.1]hept-2-yl]-5-heptenoic acid

MF: C₂₃H₂₉IO₅

FW: 512.4

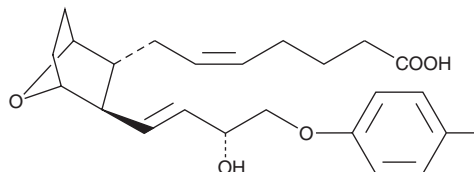
Purity: ≥95%

UV/Vis.: λ_{max}: 234 nm

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.

Laboratory Procedures

I-BOP 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 ethanol, DMSO, and dimethyl formamide purged with an inert gas can be used. The solubility of I-BOP in these solvents is approximately 100, 25, and 50 mg/ml, respectively.

I-BOP is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, the ethanolic solution of I-BOP should be diluted with the aqueous buffer of choice. The solubility of I-BOP in PBS (pH 7.2) is approximately 3 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

I-BOP is a potent agonist of TP receptor-mediated platelet aggregation and vascular smooth muscle contraction.¹ I-BOP induces platelet aggregation with EC₅₀ values of 0.34 nM and 0.174 nM at pH 7.4 and pH 6.0, respectively.²

References

1. Morinelli, T.A., Mais, D.E., Oatis, J.E., *et al.* I-BOP, the most potent radiolabelled agonist for the TXA₂/PGH₂ receptor. *Adv. Prostaglandin Thromboxane Leukotriene Res.* **20**, 102-109 (1990).
2. Mayeux, P.R., Morinelli, T.A., Williams, T.C., *et al.* Differential effect of pH on thromboxane A₂/prostaglandin H₂ receptor agonist and antagonist binding in human platelets. *J. Biol. Chem.* **266**(21), 13752-13758 (1991).

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, 06/10/2024

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