

PRODUCT INFORMATION



I-SAP

Item No. 19021

CAS Registry No.: 133538-58-6
Formal Name: (5Z)-7-[(1S,2S,3S,5R)-3-[[[4-iodophenyl)sulfonyl]amino]-6,6-dimethylbicyclo[3.1.1]hept-2-yl]-5-heptenoic acid

Synonyms: Iodophenyl sulfonyl amino pinane Thromboxane A₂, Iodophenyl sulfonyl amino pinane TXA₂

MF: C₂₂H₃₀INO₄S

FW: 531.4

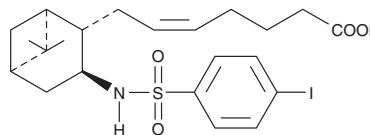
Purity: ≥98%

UV/Vis.: λ_{max}: 249 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-SAP is supplied as a solution in ethanol. To change the solvent, simply evaporate the I-SAP 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 I-SAP in these solvents is approximately 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 I-SAP is needed, it can be prepared by evaporating the ethanol and directly dissolving the neat oil in aqueous buffers. The solubility of I-SAP in PBS (pH 7.2) is approximately 100 µg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

I-SAP is a high affinity TP receptor antagonist. At physiologic pH, I-SAP produces platelet shape change, but not aggregation, with an EC₅₀ value of 9.7 nM. I-SAP binds to human platelets with the maximum binding obtained between pH 6.5 and pH 7.4. In washed human platelets, the K_d for I-SAP is 468 pM at pH 7.4 and 490 pM at pH 6.5.¹

Reference

1. Naka, M., Mais, D.E., Morinelli, T.A., *et al.* 7-[(1R,2S,3S,5R)-6,6-Dimethyl-3-(4-iodobenzenesulfonylamino)bicyclo[3.1.1]hept-2-yl]-5(Z)-heptenoic acid: A novel high-affinity radiolabeled antagonist for platelet thromboxane A₂/prostaglandin H₂ receptors. *J. Pharmacol. Exp. Ther.* **262**(2), 632-637 (1992).

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, 11/08/2021

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