

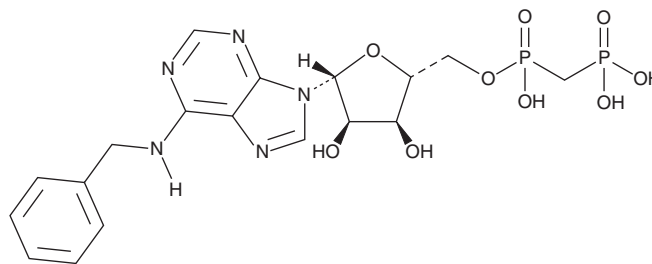
PRODUCT INFORMATION



PSB-12379

Item No. 28446

CAS Registry No.: 1802226-78-3
Formal Name: N-(phenylmethyl)-adenosine, 5'-[hydrogen P-(phosphonomethyl)phosphonate]
MF: C₁₈H₂₃N₅O₉P₂
FW: 515.4
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

PSB-12379 is supplied as a solid. A stock solution may be made by dissolving the PSB-12379 in water. The solubility of PSB-12379 in water is approximately 5 mg/ml (sonicated). We do not recommend storing the aqueous solution for more than one day.

Description

PSB-12379 is an inhibitor of ecto-5'-nucleotidase (CD73; $K_s = 2.21$ and 9.03 nM for the recombinant human and rat enzyme, respectively).¹ It is selective for CD73 over human nucleoside triphosphate diphosphohydrolase 1 (NTPDase 1), -2, and -3, ectonucleotide pyrophosphatase/phosphodiesterase family member 1 (ENPP1), -2, and -3 ($K_s = >10,000$ nM for all), as well as the purinergic P2Y₁ and P2Y₁₂ receptors at 10 μ M. PSB-12379 inhibits CD73 in isolated human serum with a K_i value of 18.8 nM.

Reference

1. Bhattarai, S., Freundlieb, M., Pippel, J., *et al.* α,β -Methylene-ADP (AOPCP) derivatives and analogues: Development of potent and selective ecto-5'-nucleotidase (CD73) inhibitors. *J. Med. Chem.* **58**, 6248-6263 (2015).

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, 12/14/2022

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