

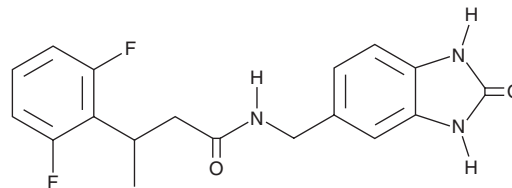
PRODUCT INFORMATION



AGPS-IN-2i

Item No. 43635

CAS Registry No.: 2316782-88-2
Formal Name: N-[(2,3-dihydro-2-oxo-1H-benzimidazol-5-yl)methyl]-2,6-difluoro- β -methyl-benzenepropanamide
Synonyms: AGPS Inhibitor 2i, alkyl-DHAP Synthase-IN-2i, alkyl-DHAP Synthase Inhibitor 2i, Alkyldihydroxyacetonephosphate Synthase-IN-21, Alkyldihydroxyacetonephosphate Synthase Inhibitor 2i



MF: C₁₈H₁₇F₂N₃O₂
FW: 345.3
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

AGPS-IN-2i is supplied as a solid. A stock solution may be made by dissolving the AGPS-IN-2i in the solvent of choice, which should be purged with an inert gas. AGPS-IN-2i is sparingly soluble (1-10 mg/ml) in DMSO.

Description

AGPS-IN-2i is an inhibitor of alkyldihydroxyacetonephosphate synthase (alkyl-DHAP synthase).¹ It decreases levels of nine ether phospholipids in 231MFP breast cancer cells when used at a concentration of 500 μ M. AGPS-IN-2i reduces migration of, and induces cytotoxicity in, 231MFP cells.

Reference

1. Stazi, G., Battistelli, C., Piano, V., *et al.* Development of alkyl glycerone phosphate synthase inhibitors: Structure-activity relationship and effects on ether lipids and epithelial-mesenchymal transition in cancer cells. *Eur. J. Med. Chem.* **163**, 722-735 (2019).

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, 09/02/2025

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