

PRODUCT INFORMATION



OPB-171775

Item No. 45601

CAS Registry No.: 2131210-15-4
Formal Name: 6-[2,3-difluoro-4-(2-hydroxy-2-methylpropoxy) phenyl]-4,5-dihydro-5-methyl-3(2H)-pyridazinone

MF: C₁₅H₁₈F₂N₂O₃

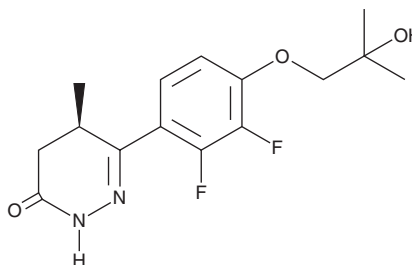
FW: 312.3

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

OPB-171775 is supplied as a solid. A stock solution may be made by dissolving the OPB-171775 in the solvent of choice, which should be purged with an inert gas. OPB-171775 is slightly soluble (0.1-1 mg/ml) in DMSO and ethanol.

Description

OPB-171775 is a molecular glue.¹ It induces the formation of a complex between phosphodiesterase 3A (PDE3A) and ribonuclease SLFN12. OBI-171775 (10 nM) induces G₀/G₁ phase cell cycle arrest, apoptosis, and necrosis in PFSK-1 primitive neuroectodermal tumor cells, with cell cycle arrest partially blocked by the general control nonderepressible 2 kinase (GCN2) inhibitor GCN2iB (Item No. 35897). It reduces tumor volume and induces tumor regression in patient-derived xenograft (PDX) mouse models of a variety of cancer types, including gastrointestinal stromal tumor (GIST), ovarian, and lung cancers. OPB-171775 (0.03 mg/kg) reduces tumor growth in a GS11353 PDX mouse model of GIST, with tumor eradication when administered in combination with imatinib (Item No. 13139).

Reference

1. Takaki, E.O., Kiyono, K., Obuchi, Y., *et al.* A PDE3A-SLFN12 molecular glue exhibits significant antitumor activity in TKI-resistant gastrointestinal stromal tumors. *Clin. Cancer Res.* **30(16)**, 3603-3621 (2024).

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, 05/04/2026

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