

PRODUCT INFORMATION

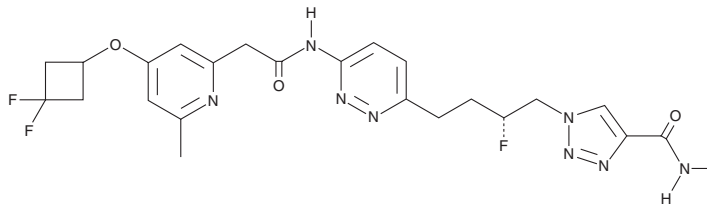


IPN60090

Item No. 35855

CAS Registry No.: 1853164-83-6
Formal Name: 4-[(3,3-difluorocyclobutyl)oxy]-N-[6-[(3R)-3-fluoro-4-[4-[(methylamino)carbonyl]-1H-1,2,3-triazol-1-yl]butyl]-3-pyridazinyl]-6-methyl-2-pyridineacetamide

Synonym: IACS-6274
MF: C₂₄H₂₇F₃N₈O₃
FW: 532.5
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

IPN60090 is supplied as a solid. A stock solution may be made by dissolving the IPN60090 in the solvent of choice, which should be purged with an inert gas. IPN60090 is soluble in methanol and DMSO.

Description

IPN60090 is an inhibitor of glutaminase 1 (GLS1; IC₅₀ = 0.031 μM for the glutaminase C isoform).¹ It is selective for GLS1 over GLS2 (IC₅₀ = >50 μM), as well as panels of 80 ion channels and receptors and 97 kinases at 10 μM. IPN60090 also binds to zinc finger SWIM domain-containing protein 4 (ZSWIM4; K_d = 2.98 μM).² It decreases the viability of A549 non-small cell lung cancer (NSCLC) cells (IC₅₀ = 26 nM) and decreases glycine levels in OVCAR-8 and SKOV3 ovarian cancer cells.^{1,2} IPN60090 (30 mg/kg, p.o.), in combination with the DNA-crosslinking agent carboplatin (Item No. 13112), decreases tumor weight and volume in a SKOV3 mouse xenograft model.²

References

1. Soth, M.J., Le, K., Di Francesco, M.E., *et al.* Discovery of IPN60090, a clinical stage selective glutaminase-1 (GLS-1) inhibitor with excellent pharmacokinetic and physicochemical properties. *J. Med. Chem.* **63(21)**, 12957-12977 (2020).
2. Gong, K., Huang, Y., Zheng, Y., *et al.* ZSWIM4 inhibition improves chemosensitivity in epithelial ovarian cancer cells by suppressing intracellular glycine biosynthesis. *J. Transl. Med.* **22(1)**, 192 (2021).

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, 03/26/2024

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