

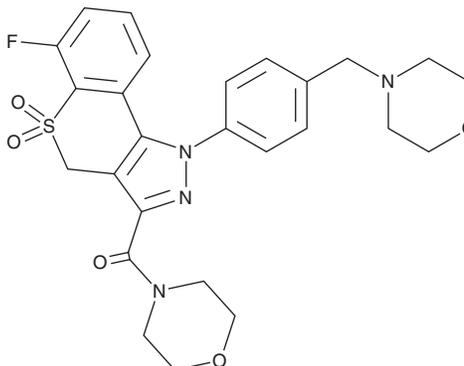
PRODUCT INFORMATION



Roginolisib

Item No. 44445

CAS Registry No.: 1305267-37-1
Formal Name: [6-fluoro-1,4-dihydro-1-[4-(4-morpholinylmethyl)phenyl]-5,5-dioxido[1]benzothiopyrano[4,3-c]pyrazol-3-yl]-4-morpholinyl-methanone
Synonyms: IOA-244, MSC2360844
MF: C₂₆H₂₇FN₄O₅S
FW: 526.6
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

Roginolisib is supplied as a solid. A stock solution may be made by dissolving the roginolisib in the solvent of choice, which should be purged with an inert gas. Roginolisib is slightly soluble (0.1-1 mg/ml) in acetonitrile:water (1:1).

Description

Roginolisib is an inhibitor of PI3K δ (IC₅₀ = 0.019 μ M).¹ It is selective for PI3K δ over PI3K α , PI3K β , PI3K γ , VPS34, and PI3K ζ (IC₅₀s = 0.43, 1.5, 10.3, 5.6, and 5.9 μ M, respectively). Roginolisib inhibits the growth of SU-DHL-6 B cell lymphoma cells that endogenously express PI3K δ (GI₅₀ = 0.24 μ M) but not cancer cell lines endogenously expressing PI3K α , PI3K β , or PI3K γ . *In vivo*, roginolisib (30 mg/kg twice daily) alone or in combination with an anti-PD-1 antibody slows tumor growth and decreases tumor volume in a murine Lewis lung carcinoma model.

Reference

1. Johnson, Z., Tarantelli, C., Civanelli, E., et al. IOA-244 is a non-ATP-competitive, highly selective, tolerable PI3K delta inhibitor that targets solid tumors and breaks immune tolerance. *Cancer Res. Commun.* **3(4)**, 576-591 (2023).

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, 12/22/2025

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