

PRODUCT INFORMATION

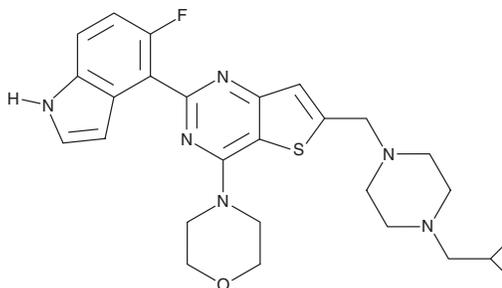


PI-3065

Item No. 9002394

CAS Registry No.: 955977-50-1
Formal Name: 6-[[4-(cyclopropylmethyl)-1-piperazinyl]methyl]-2-(5-fluoro-1H-indol-4-yl)-4-(4-morpholinyl)-thieno[3,2-d]pyrimidine

MF: C₂₇H₃₁FN₆OS
FW: 506.6
Purity: ≥98%
UV/Vis.: λ_{max}: 209, 299 nm
Supplied as: A crystalline 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

PI-3065 is supplied as a crystalline solid. A stock solution may be made by dissolving the PI-3065 in the solvent of choice. PI-3065 is soluble in the organic solvent DMSO, which should be purged with an inert gas, at a concentration of approximately 50 mg/ml. Slight warming of the DMSO may be necessary.

Description

PI-3065 is an inhibitor of PI3K p110δ (IC₅₀ = 15 nM).¹ It is greater than 100-fold selective for p110δ over p110α, p110β, p110γ, DNA protein kinase (DNA-PK), and mTOR. PI-3065 (75 mg/kg) decreases tumor weight and inhibits metastasis in a 4T1 murine mammary carcinoma model.

The information regarding the kinase IC₅₀ values and xenograft model was drawn from a paper that has been retracted; however, the information specified in the retraction statement has not been included.¹

Reference

1. Ali, K., Soond, D.R., Pineiro, R., *et al.* Inactivation of PI3K p110δ breaks regulatory T-cell-mediated immune tolerance to cancer. *Nature* **510(7505)**, 407-411 (2014).

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, 10/24/2023

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