

PRODUCT INFORMATION



PI3-Kinase Inhibitor

Item No. 43486

CAS Registry No.: 2504036-13-7
Formal Name: N-[5-[2-[(1S)-1-cyclopropylethyl]-2,3-dihydro-4-(4-morpholinyl)-3-oxo-1H-pyrrolo[3,4-c]pyridin-6-yl]-4-methyl-2-thiazolyl]-acetamide

Synonyms: PI3K Inhibitor, Phosphatidylinositol 3-Kinase Inhibitor

MF: C₂₂H₂₇N₅O₃S

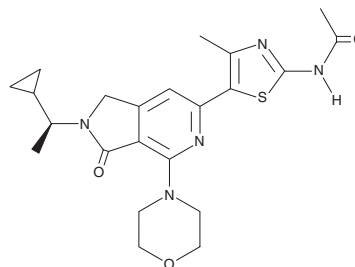
FW: 441.6

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

PI3-Kinase inhibitor is supplied as a solid. A stock solution may be made by dissolving the PI3-kinase inhibitor in the solvent of choice, which should be purged with an inert gas. PI3-Kinase inhibitor is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

Description

PI3-Kinase inhibitor is an inhibitor of PI3K γ (IC₅₀ = 1.5 nM).¹ It is selective for PI3K γ over PI3K α , PI3K β , and PI3K δ (IC₅₀s = 3,333, >10,000, and >1,000 nM, respectively). PI3-Kinase inhibitor (100 mg/kg) inhibits tumor growth in a 4T1 murine mammary cancer model. It decreases disease severity in a mouse model of experimental autoimmune encephalomyelitis (EAE) when administered at doses of 20 or 80 mg/kg.

Reference

1. Hao, X. Phosphatidylinositol 3-kinase inhibitors. **WO2020/210379A1** (2020).

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, 07/09/2025

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