

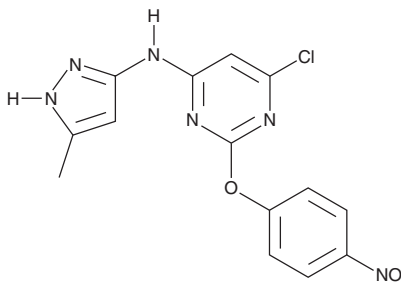
PRODUCT INFORMATION



Autophinib

Item No. 27320

CAS Registry No.: 1644443-47-9
Formal Name: 6-chloro-N-(5-methyl-1H-pyrazol-3-yl)-2-(4-nitrophenoxy)-4-pyrimidinamine
MF: C₁₄H₁₁ClN₆O₃
FW: 346.7
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

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

Description

Autophinib is an ATP-competitive inhibitor of the PI3K VPS34 ($IC_{50} = 19 \text{ nM}$).¹ It inhibits autophagy induced by the mTOR inhibitor rapamycin (Item No. 13346) or by amino acid starvation (IC_{50} s = 40 and 90 nM, respectively) in MCF-7 cells. Autophinib also induces apoptosis in amino acid-starved MCF-7 cells with an EC_{50} value of 234 nM.

Reference

1. Robke, L., Laraia, L., Carnero Corrales, M.A., *et al.* Phenotypic identification of a novel autophagy inhibitor chemotype targeting lipid kinase VPS34. *Angew. Chem. Int. Ed. Engl.* **56(28)**, 8153-8157 (2017).

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, 12/16/2022

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