

PRODUCT INFORMATION

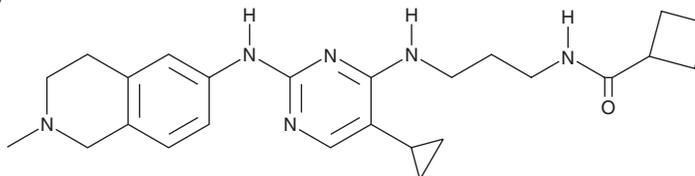


MRT68921

Item No. 19905

CAS Registry No.: 1190379-70-4
Formal Name: N-[3-[[5-cyclopropyl-2-[(1,2,3,4-tetrahydro-2-methyl-6-isoquinoliny)amino]-4-pyrimidinyl]amino]propyl]-cyclobutanecarboxamide

MF: C₂₅H₃₄N₆O
FW: 434.6
Purity: ≥98%
UV/Vis.: λ_{max}: 208, 262 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

MRT68921 is supplied as a crystalline solid. A stock solution may be made by dissolving the MRT68921 in the solvent of choice, which should be purged with an inert gas. MRT68921 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of MRT68921 in these solvents is approximately 2, 3, and 5 mg/ml, respectively.

MRT68921 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, MRT68921 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. MRT68921 has a solubility of approximately 0.16 mg/ml in a 1:5 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

MRT68921 is an inhibitor of ULK1 and ULK2 (IC₅₀s = 2.9 and 1.1 nM, respectively).¹ Through its effects on ULK1, MRT68921 blocks autophagy in cells, driving the accumulation of stalled early autophagosomal structures.¹

Reference

1. Petherick, K.J., Conway, O.J.L., Mpamhanga, C., *et al.* Pharmacological inhibition of ULK1 kinase blocks mammalian target of rapamycin (mTOR)-dependent autophagy. *J. Biol. Chem.* **290**(18), 11376-11383 (2015).

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, 09/30/2022

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