

PRODUCT INFORMATION

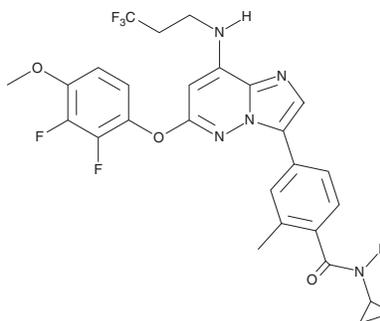


BAY 12-17389

Item No. 21660

CAS Registry No.: 1554458-53-5
Formal Name: N-cyclopropyl-4-[6-(2,3-difluoro-4-methoxyphenoxy)-8-[(3,3,3-trifluoropropyl)amino]imidazo[1,2-b]pyridazin-3-yl]-2-methyl-benzamide

MF: C₂₇H₂₄F₅N₅O₃
FW: 561.5
Purity: ≥95%
UV/Vis.: λ_{max}: 271, 304 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

BAY 12-17389 is supplied as a crystalline solid. A stock solution may be made by dissolving the BAY 12-17389 in the solvent of choice, which should be purged with an inert gas. BAY 1217389 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of BAY 12-17389 in these solvents is approximately 15 and 25 mg/ml, respectively.

BAY 12-17389 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, BAY 12-17389 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. BAY 12-17389 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

BAY 12-17389 is an inhibitor of monopolar spindle 1 (Mps1; IC₅₀ = 0.63 nM).¹ It is selective for Mps1 over a panel of kinases at 10 nM. BAY 12-17389 promotes cell-cycle progression into mitosis in metaphase-arrested HeLa cells (IC₅₀ = 11 nM). BAY 12-17389 (1.5 mg/kg) enhances paclitaxel-induced inhibition of tumor growth in a SUM149 triple-negative breast cancer (TNBC) mouse xenograft model.

Reference

1. Wengner, A.M., Siemeister, G., Koppitz, M., *et al.* Novel Mps1 kinase inhibitors with potent antitumor activity. *Mol. Cancer Ther.* **15**(4), 583-592 (2016).

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, 01/20/2023

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