

PRODUCT INFORMATION

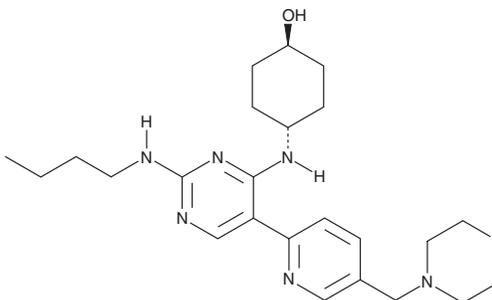


UNC2250

Item No. 17703

CAS Registry No.: 1493694-70-4
Formal Name: *trans*-4-[[2-(butylamino)-5-[5-(4-morpholinylmethyl)-2-pyridinyl]-4-pyrimidinyl]amino]-cyclohexanol

MF: C₂₄H₃₆N₆O₂
FW: 440.6
Purity: ≥98%
UV/Vis.: λ_{max}: 232, 274, 332 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.

Description

Tyros3, Axl, and Mer are members of the TAM family of receptor tyrosine kinases that have signaling roles in normal and malignant cells.¹ UNC2250 is an inhibitor of Mer kinase activity, blocking the steady-state phosphorylation of endogenous Mer in 697 B-ALL cells with an IC₅₀ value of 9.8 nM.² It is selective for Mer over a panel of both tyrosine and serine-threonine kinases, including Tyros3 and Axl. UNC2250 also inhibits ligand-dependent phosphorylation of a chimeric EGFR-Mer protein and blocks colony formation in soft agar cultures of BT-12 rhabdoid tumor and Colo699 NSCLC cell lines.² UNC2250 displays good pharmacokinetic properties in mice following intravenous or oral administration.²

Reference

1. Verma, A., Warner, S.L., Vankayalapati, H., *et al.* Targeting Axl and Mer kinases in cancer. *Mol. Cancer Ther.* **10**(10), 1763-1773 (2016).
2. Zhang, W., Zhang, D., Stashko, M.A., *et al.* Pseudo-cyclization through intramolecular hydrogen bond enables discovery of pyridine substituted pyrimidines as new Mer kinase inhibitors. *J. Med. Chem.* **56**(23), 9683-9692 (2013).

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, 12/06/2022

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