

PRODUCT INFORMATION

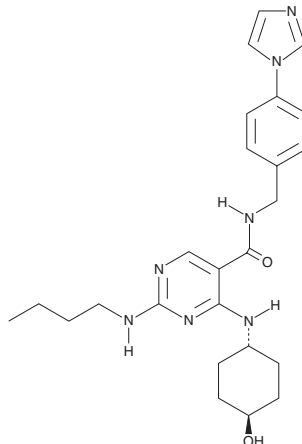


UNC2881

Item No. 29012

CAS Registry No.: 1493764-08-1
Formal Name: 2-(butylamino)-4-[(*trans*-4-hydroxycyclohexyl)amino]-N-[[4-(1H-imidazol-1-yl)phenyl]methyl]-5-pyrimidinecarboxamide

MF: C₂₅H₃₃N₇O₂
FW: 463.6
Purity: ≥98%
UV/Vis.: λ_{max}: 232, 263, 309 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

UNC2881 is supplied as a crystalline solid. A stock solution may be made by dissolving the UNC2881 in the solvent of choice, which should be purged with an inert gas. UNC2881 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of UNC2881 in ethanol is approximately 1 mg/ml and approximately 30 mg/ml in DMSO and DMF.

UNC2881 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, UNC2881 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. UNC2881 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

UNC2881 is an inhibitor of Mer (IC₅₀ = 4.3 nM), a member of the TAM family of receptor tyrosine kinases.¹ It is selective for Mer over the remaining TAM family members Axl and TYRO3 (IC₅₀s = 360 and 250 nM, respectively). UNC2881 inhibits phosphorylation of Mer in 697 B-ALL acute lymphoblastic leukemia cells (IC₅₀ = 21.9 nM). It inhibits platelet aggregation and ATP release induced by type I equine fibrillar collagen in isolated human platelet-rich plasma when used at a concentration of 3 μM.

Reference

1. Zhang, W., McIver, A.L., Stashko, M.A., *et al.* Discovery of Mer specific tyrosine kinase inhibitors for the treatment and prevention of thrombosis. *J. Med. Chem.* **56**(23), 9693-9700 (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, 11/02/2022

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