

PRODUCT INFORMATION

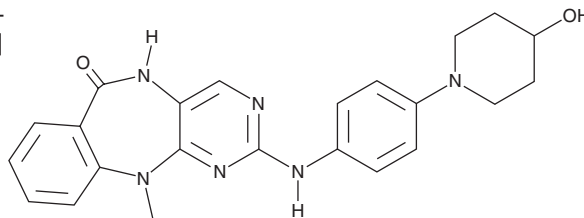


XMD16-5

Item No. 23353

CAS Registry No.: 1345098-78-3
Formal Name: 5,11-dihydro-2-[[4-(4-hydroxy-1-piperidiny)phenyl]amino]-11-methyl-6H-pyrimido[4,5-b][1,4]benzodiazepin-6-one

MF: C₂₃H₂₄N₆O₂
FW: 416.5
Purity: ≥98%
UV/Vis.: λ_{max}: 297 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

XMD16-5 is supplied as a crystalline solid. A stock solution may be made by dissolving the XMD16-5 in the solvent of choice, which should be purged with an inert gas. XMD16-5 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of XMD16-5 in these solvents is approximately 33 mg/ml.

XMD16-5 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, XMD16-5 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. XMD16-5 has a solubility of approximately 0.125 mg/ml in a 1:7 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

XMD16-5 is an inhibitor of tyrosine kinase non-receptor 2 (TNK2; IC₅₀ = 0.38 μM), which contains mutations in certain cancer cells.¹ It also decreases viability of Ba/F3 cells expressing the oncogenic TNK2 D163E and R806Q mutant proteins (IC₅₀s = 16 and 77 nM, respectively) without affecting viability of parental cell lines up to a concentration of approximately 1 μM.

Reference

1. Maxson, J.E., Abel, M.L., Wang, J., *et al.* Identification and characterization of tyrosine kinase nonreceptor 2 mutations in leukemia through integration of kinase inhibitor screening and genomic analysis. *Cancer Res.* **76**(1), 127-138 (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, 11/30/2022

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