

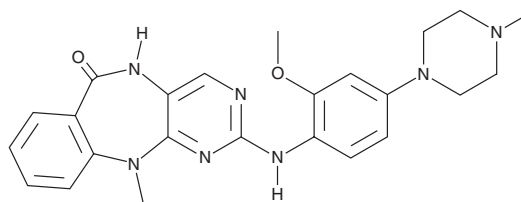
PRODUCT INFORMATION



XMD8-87

Item No. 23354

CAS Registry No.: 1234480-46-6
Formal Name: 5,11-dihydro-2-[[2-methoxy-4-(4-methyl-1-piperazinyl)phenyl]amino]-11-methyl-6H-pyrimido[4,5-b][1,4]benzodiazepin-6-one
MF: C₂₄H₂₇N₇O₂
FW: 445.5
Purity: ≥98%
UV/Vis.: λ_{max}: 204, 295 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

XMD8-87 is supplied as a crystalline solid. A stock solution may be made by dissolving the XMD8-87 in the solvent of choice, which should be purged with an inert gas. XMD8-87 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of XMD8-87 in these solvents is approximately 1 mg/ml.

Description

XMD8-87 is an inhibitor of tyrosine kinase non-receptor 2 (TNK2) with an IC₅₀ value of 1.9 μM in HEK293T cells expressing human TNK2.¹ It is selective for TNK2 over a panel of approximately 100 kinases with >90 and <50% inhibition for TNK2 and all other kinases tested, respectively. XMD8-87 inhibits growth of Ba/F3 cells expressing constitutively active mutant TNK2s (IC₅₀s = 0.1-77 nM) commonly found in leukemias but has no effect on cells expressing wild-type TNK2 (IC₅₀ = >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, 12/16/2022

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