

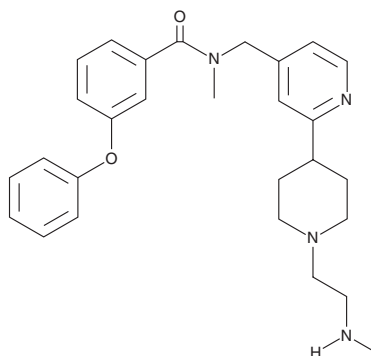
PRODUCT INFORMATION



TP-064

Item No. 20256

CAS Registry No.: 2080306-20-1
Formal Name: N-methyl-N-[[2-[1-[2-(methylamino)ethyl]-4-piperidinyl]-4-pyridinyl]methyl]-3-phenoxy-benzamide
MF: C₂₈H₃₄N₄O₂
FW: 458.6
Purity: ≥98%
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

TP-064 is supplied as a crystalline solid. A stock solution may be made by dissolving the TP-064 in the solvent of choice, which should be purged with an inert gas. TP-064 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of TP-064 in these solvents is approximately 30 mg/ml.

Description

TP-064 is a potent inhibitor of protein arginine methyltransferase 4 (PRMT4; IC₅₀ < 10 nM) that is greater than 100-fold selective over other histone methyltransferases and non-epigenetic targets. It also inhibits methylation of the RNA polymerase II mediator complex subunit 12 (MED12; IC₅₀ = 43 nM) *in vitro*. For more information on TP-064 please visit the Structural Genomics Consortium (SGC). The negative control, TP-064N, for TP-064 is also available exclusively through the SGC.

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, 11/11/2022

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