

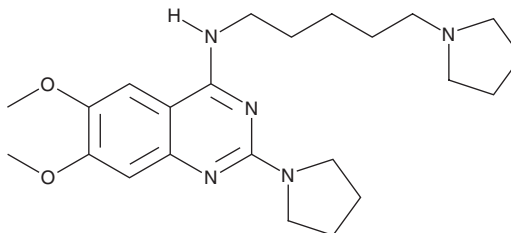
PRODUCT INFORMATION



UNC0379

Item No. 16400

CAS Registry No.: 1620401-82-2
Formal Name: 6,7-dimethoxy-2-(1-pyrrolidinyl)-
N-[5-(1-pyrrolidinyl)pentyl]-4-
quinazolinamine
MF: C₂₃H₃₅N₅O₂
FW: 413.6
Purity: ≥95%
UV/Vis.: λ_{max}: 210, 249, 345 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

UNC0379 is supplied as a crystalline solid. A stock solution may be made by dissolving the UNC0379 in the solvent of choice. UNC0379 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of UNC0379 in ethanol is approximately 10 mg/ml and approximately 20 mg/ml in DMSO and DMF.

UNC0379 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, UNC0379 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. UNC0379 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

SET domain-containing protein 8 (SET8) is a methyltransferase that selectively mono-methylates histone H4 at lysine residue 20, an event proven to have an important role in chromatin structure and transcriptional activation. It is also a regulator of p53, mono-methylating lysine 382 of the tumor suppressor. UNC0379 is a substrate-competitive inhibitor of the lysine methyltransferase SET8 (IC₅₀ = 7.3 μM; K_d = 18.3 μM).¹ It is selective for SET8 over 15 other methyltransferases, including G9a and GLP (IC₅₀s = >100 μM).¹

Reference

1. Ma, A., Yu, W., Li, F., *et al.* Discovery of a selective, substrate-competitive inhibitor of the lysine methyltransferase SETD8. *J. Med. Chem.* (2014).

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, 11/16/2022

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