

PRODUCT INFORMATION

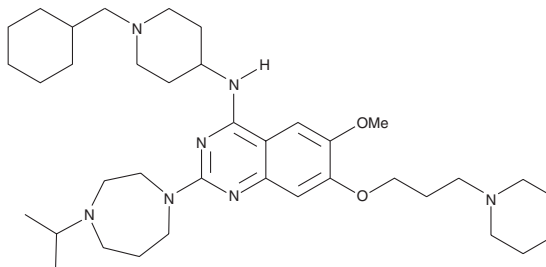


UNC0631

Item No. 11084

CAS Registry No.: 1320288-19-4
Formal Name: N-[1-(cyclohexylmethyl)-4-piperidyl]-2-[hexahydro-4-(1-methylethyl)-1H-1,4-diazepin-1-yl]-6-methoxy-7-[3-(1-piperidyl)propoxy] 4-quinazolinamine

MF: C₃₇H₆₁N₇O₂
FW: 635.9
Purity: ≥98%
UV/Vis.: λ_{max}: 251, 348 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

UNC0631 is supplied as a crystalline solid. A stock solution may be made by dissolving the UNC0631 in the solvent of choice, which should be purged with an inert gas. UNC0631 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of UNC0631 in DMSO is approximately 5 mg/ml and approximately 30 mg/ml in ethanol and DMF.

UNC0631 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, UNC0631 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. UNC0631 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

UNC0631 is a potent and selective inhibitor of G9a activity *in vitro* (IC₅₀ = 4 nM) and G9a/GLP-mediated dimethylation of histone 3 on lysine 9 in MDA-MB-231 cells (IC₅₀ = 25 nM).¹ It potently inhibits G9a/GLP activity in a variety of cancer cell lines as well as in the human fetal lung IMR-90 cell line.¹

Reference

1. Liu, F., Barysyt-Lovejoy, D., Allali-Hassani, A., *et al.* Optimization of cellular activity of G9a inhibitors 7-aminoalkoxy-quinazolines. *J. Med. Chem.* **54**(17), 6139-6150 (2011).

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, 11/14/2022

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