
UNC0638 hydrate

(Catalog No. M63000)

Alternative Names

2-cyclohexyl-6-methoxy-N-[1-(1-methylethyl)-4-piperidinyl]-7-[3-(1-pyrrolidinyl)propoxy]-4-quinazolinamine

Description

1,2,3 UNC0638 is a potent G9a HMTase inhibitor, exhibiting an IC₅₀ value of <15 nM in vitro. UNC0638 also inhibits GLP, a closely-related H3K9 HMTase, with an IC₅₀ value of 19 nM, but is more than 10,000-fold selective against SET7/9 (a H3K4 HMTase), SET8 (a H4K20 HMTase), PRMT3, and SUV39H2.

CAS Number

1255580-76-7

Chemical Formula

C₃₀H₄₇N₅O₂

Molecular Weight

509.7

Size

1 mg

Purity

≥98%

Solubility

Soluble in DMSO

Storage Temp

-20°C

Ordering Information:

Products	Size	Cat. No.
UNC0638 hydrate	1 mg	M63000 -1

This product is for research purposes only. Not intended for use in diagnostic procedures.