

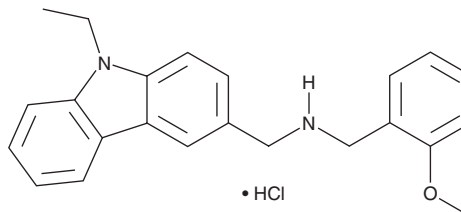
PRODUCT INFORMATION



HLCL-61 (hydrochloride)

Item No. 19897

CAS Registry No.: 1158279-20-9
Formal Name: 9-ethyl-N-[(2-methoxyphenyl)methyl]-9H-carbazole-3-methanamine, monohydrochloride
MF: $C_{23}H_{24}N_2O \cdot HCl$
FW: 380.9
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 239, 267, 297 nm
Supplied as: A crystalline solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

HLCL-61 (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the HLCL-61 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. HLCL-61 (hydrochloride) is soluble in organic solvents such as ethanol and DMSO. The solubility of HLCL-61 (hydrochloride) in these solvents is approximately 1 mg/ml.

Description

HLCL-61 is an inhibitor of protein arginine methyltransferase 5 (PRMT5) that inhibits the growth of multiple acute myeloid leukemia (AML) cell lines and patient-derived tumor samples ($IC_{50}s = 7.21$ - 21.46 and 3.98 - $8.72 \mu M$, respectively).¹ It is selective for PRMT5, lacking activity against PRMT1, PRMT4, and PRMT7 in an enzyme assay. HLCL-61 induces myeloid differentiation of THP-1 cells and increases CD11b expression in a dose-dependent manner. It also increases expression of miR-29b mRNA resulting in a 4-fold decrease in FLT3 activity in THP-1 cells expressing an FLT3 luciferase reporter.

Reference

1. Tarighat, S.S., Santhanam, R., Frankhouser, D., *et al.* The dual epigenetic role of PRMT5 in acute myeloid leukemia: Gene activation and repression via histone arginine methylation. *Leukemia* **30**(4), 789-799 (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, 10/26/2022

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