

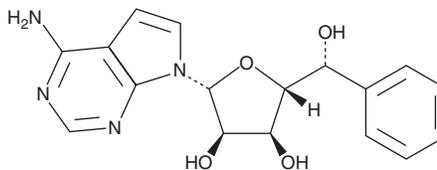
PRODUCT INFORMATION



LLY-283

Item No. 21596

CAS Registry No.: 2040291-27-6
Formal Name: (7-[(5R)-5-C-phenyl-β-D-ribofuranosyl]-7H-pyrrolo[2,3-d]pyrimidin-4-amine
MF: C₁₇H₁₈N₄O₄
FW: 342.4
Purity: ≥98%
UV/Vis.: λ_{max}: 230, 274 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

LLY-283 is supplied as a crystalline solid. A stock solution may be made by dissolving the LLY-283 in the solvent of choice. LLY-283 is soluble in organic solvents such as ethanol and DMSO, which should be purged with an inert gas. The solubility of LLY-283 in these solvents is approximately 60 mg/ml.

Description

LLY-283 is an inhibitor of protein arginine methyltransferase 5 (PRMT5; IC₅₀ = 22 nM).¹ It is selective for PRMT5 over a panel of methyltransferases, including PRMT4, -6, and -7 at 1 μM. It reduces symmetric demethylation of SmBB' in MCF-7 cells (IC₅₀ = 25 nM) and inhibits proliferation of various breast, gastric, hematological, lung, skin, and ovarian cancer cell lines (IC₅₀s = 3-30 nM). LLY-283 (20 mg/kg) inhibits tumor growth in an A375 mouse xenograft model. It also disrupts alternative splicing events in, and inhibits proliferation and self-renewal of, patient-derived glioblastoma stem cells and increases survival in an orthotopic patient-derived xenograft (PDX) mouse model.² See the Structural Genomics Consortium (SGC) website for more information.

References

1. Bonday, Z.Q., Cortez, G.S., Grogan, M.J., *et al.* LLY-283, a potent and selective inhibitor of arginine methyltransferase 5, PRMT5, with antitumor activity. *ACS Med. Chem. Lett.* **9(7)**, 612-617 (2018).
2. Sachamitr, P., Ho, J.C., Ciamponi, F.E., *et al.* PRMT5 inhibition disrupts splicing and stemness in glioblastoma. *Nat Commun.* **12(1)**, 979 (2021).

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, 11/10/2022

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