

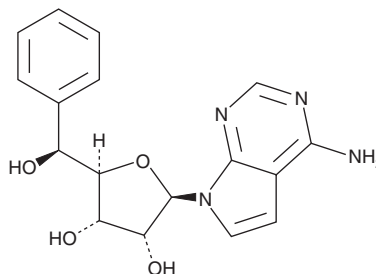
PRODUCT INFORMATION



LLY-284

Item No. 43141

CAS Registry No.: 2226515-75-7
Formal Name: 7-[(5S)-5-C-phenyl-β-D-ribofuranosyl]-7H-pyrrolo[2,3-d]pyrimidin-4-amine
MF: C₁₇H₁₈N₄O₄
FW: 342.4
Purity: ≥98%
Supplied as: A 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-284 is supplied as a solid. A stock solution may be made by dissolving the LLY-284 in the solvent of choice, which should be purged with an inert gas. LLY-284 is soluble (≥10 mg/ml) in ethanol and DMSO.

Description

LLY-284 is a diastereomer of the protein arginine methyltransferase 5 (PRMT5) inhibitor LLY-283 (Item No. 21596).¹ It weakly inhibits PRMT5 (IC₅₀ = 1,074 nM). LLY-284 (100 μM) inhibits the asexual growth of *P. falciparum* *in vitro*.²

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. Min, H., Lucky, A.B., Madsen, J.J., *et al.* Onametostat, a PfPRMT5 inhibitor, exhibits antimalarial activity to *Plasmodium falciparum*. *Antimicrob. Agents Chemother.* **68**(10), e0017624 (2024).

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, 04/17/2025

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