

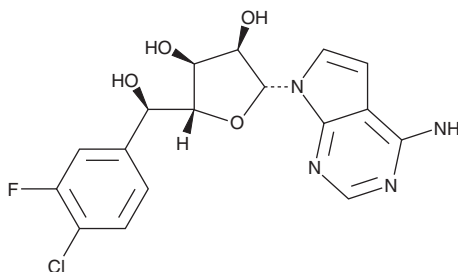
PRODUCT INFORMATION



PF-06855800

Item No. 43630

CAS Registry No.: 1989620-04-3
Formal Name: 7-[(5R)-5-C-(4-chloro-3-fluorophenyl)- β -D-ribofuranosyl]-7H-pyrrolo[2,3-d]pyrimidin-4-amine
MF: C₁₇H₁₆ClFN₄O₄
FW: 394.8
Purity: \geq 98%
Supplied as: A solid
Storage: -20°C
Stability: \geq 4 years



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

Laboratory Procedures

PF-06855800 is supplied as a solid. A stock solution may be made by dissolving the PF-06855800 in the solvent of choice, which should be purged with an inert gas. PF-06855800 is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

Description

PF-06855800 is an inhibitor of protein arginine methyltransferase 5 (PRMT5; $K_i = 11$ pM).¹ It is selective for PRMT5 over a panel of 22 additional protein methyltransferases and a panel of 40 kinases at 10 μ M. PF-06855800 inhibits the proliferation of A2780 ovarian cancer cells ($IC_{50} = 3.3$ nM). It also inhibits the asexual growth of *P. falciparum* *in vitro* when used at a concentration of 100 μ M.²

References

- Jensen-Pergakes, K., Tatlock, J., Maegley, K.A., *et al.* SAM-competitive PRMT5 inhibitor PF-06939999 demonstrates antitumor activity in splicing dysregulated NSCLC with decreased liability of drug resistance. *Mol. Cancer Ther.* **21(1)**, 3-15 (2022).
- 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, 07/24/2025

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