

# PRODUCT INFORMATION



## PF-06939999 (hydrochloride)

Item No. 44357

**Formal Name:** (1S,2S,3S,5R)-3-[[6-(difluoromethyl)-5-fluoro-1,2,3,4-tetrahydro-8-isoquinolinyl]oxy]-5-(4-methyl-7H-pyrrolo[2,3-d]pyrimidin-7-yl)-1,2-cyclopentanediol, dihydrochloride

**Synonyms:** PRMT5-IN-3, PRMT5 Inhibitor 3, Protein Arginine Methyltransferase 5-IN-3, Protein Arginine Methyltransferase 5 Inhibitor 3

**MF:** C<sub>22</sub>H<sub>23</sub>F<sub>3</sub>N<sub>4</sub>O<sub>3</sub> • 2HCl

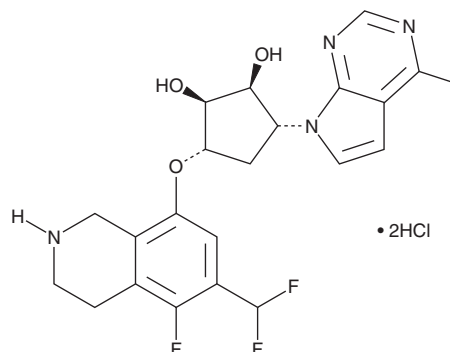
**FW:** 521.3

**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

PF-06939999 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the PF-06939999 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. PF-06939999 (hydrochloride) is slightly soluble (0.1-1 mg/ml) in acetonitrile. It is also slightly soluble (0.1-1 mg/ml) in water. We do not recommend storing the aqueous solution for more than one day.

### Description

PF-06939999 is an inhibitor of protein arginine methyltransferase 5 (PRMT5;  $K_i = <5 \text{ pM}$ ).<sup>1</sup> It is selective for PRMT5 over a panel of 22 additional protein methyltransferases and a panel of 40 kinases at 10  $\mu\text{M}$ . PF-06939999 decreases global symmetric dimethyl arginine levels ( $\text{IC}_{50} = 1.1 \text{ nM}$ ) but does not affect histone methylation levels in A-427 non-small cell lung cancer (NSCLC) cells. *In vivo*, PF-06939999 (3, 10, and 30 mg/kg) reduces tumor volume in an A-427 mouse xenograft model.

### Reference

1. 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).

**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, 09/17/2025

### 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