

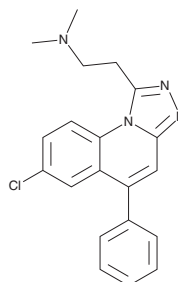
PRODUCT INFORMATION



PF-9366

Item No. 30782

CAS Registry No.: 72882-78-1
Formal Name: 7-chloro-N,N-dimethyl-5-phenyl-[1,2,4]triazolo[4,3-a]quinoline-1-ethanamine
MF: C₂₀H₁₉ClN₄
FW: 350.8
Purity: ≥98%
UV/Vis.: λ_{max}: 233 nm
Supplied as: A solid
Storage: -20°C
Stability: ≥2 years



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

Laboratory Procedures

PF-9366 is supplied as a solid. A stock solution may be made by dissolving the PF-9366 in the solvent of choice, which should be purged with an inert gas. PF-9366 is soluble in organic solvents such as ethanol and dimethyl formamide. The solubility of PF-9366 in these solvents is approximately 2 and 0.5 mg/ml, respectively.

PF-9366 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, PF-9366 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. PF-9366 has a solubility of approximately 0.2 mg/ml in a 1:4 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

PF-9366 is an inhibitor of methionine adenosyltransferase 2A (Mat2A; IC₅₀ = 420 nM for S-adenosyl-L-methionine (SAM) production in an enzyme assay).¹ It is selective for Mat2A over a panel of neurotransmitter transporters, phosphodiesterases, ion channels, and G protein-coupled receptors, as well as a panel of 39 kinases at 10 μM. PF-9366 inhibits SAM production in H520 lung carcinoma cells (IC₅₀ = 1.2 μM) and Huh7 hepatocellular carcinoma cells (IC₅₀ = 225 nM).

Reference

1. Quinlan, C.L., Kaiser, S.E., Bolaños, B., *et al.* Targeting S-adenosylmethionine biosynthesis with a novel allosteric inhibitor of Mat2A. *Nat. Chem. Biol.* **13**(7), 785-792 (2017).

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, 05/28/2020

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