

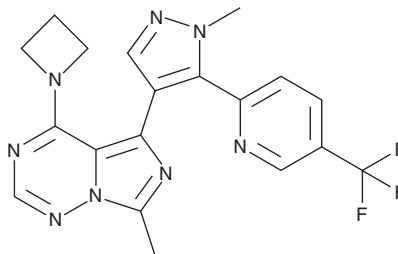
PRODUCT INFORMATION



PF-05180999

Item No. 30374

CAS Registry No.: 1394033-54-5
Formal Name: 4-(1-azetidinyl)-7-methyl-5-[1-methyl-5-[5-(trifluoromethyl)-2-pyridinyl]-1H-pyrazol-4-yl]-imidazo[5,1-f][1,2,4]triazine
MF: C₁₉H₁₇F₃N₈
FW: 414.4
Purity: ≥98%
UV/Vis.: λ_{max}: 258 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

PF-05180999 is supplied as a crystalline solid. A stock solution may be made by dissolving the PF-05180999 in the solvent of choice, which should be purged with an inert gas. PF-05180999 is soluble in the organic solvent DMSO at a concentration of approximately 10 mM.

Description

PF-05180999 is a phosphodiesterase 2A (PDE2A) inhibitor (IC₅₀ = 1.6 nM).¹ It is greater than 2000-fold selective for PDE2A over a panel of ten additional PDEs. PF-05180999 increases striatal, cortical, and hippocampal cGMP levels in mice in a dose-dependent manner. *In vivo*, it inhibits disruption of low frequency cortical δ oscillation induced by the NMDA antagonist MK-801 in rat brain when administered at a dose of 0.03 mg/kg. PF-05180999 (0.1 and 0.32 mg/kg) inhibits ketamine-induced increases in the number of working memory errors in a radial arm maze task in rats.

Reference

1. Helal, C.J., Arnold, E., Boyden, T., *et al.* Identification of a potent, highly selective, and brain penetrant phosphodiesterase 2A inhibitor clinical candidate. *J. Med. Chem.* **61**(3), 1001-1018 (2018).

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, 11/08/2022

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