

PRODUCT INFORMATION



PF-00489791

Item No. 40464

CAS Registry No.: 853003-48-2
Formal Name: 1-(2-ethoxyethyl)-5-(ethylmethylamino)-7-[(4-methyl-2-pyridinyl)amino]-N-(methylsulfonyl)-1H-pyrazolo[4,3-d]pyrimidine-3-carboxamide

Synonyms: PF-489791, UK 489791

MF: C₂₀H₂₈N₈O₄S

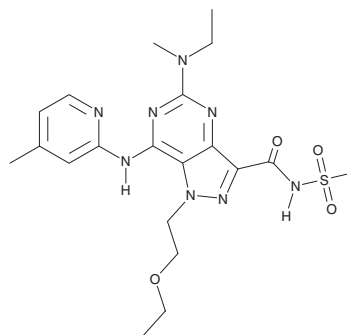
FW: 476.6

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-00489791 is supplied as a solid. A stock solution may be made by dissolving the PF-00489791 in the solvent of choice, which should be purged with an inert gas. PF-00489791 is soluble in acetonitrile. PF-00489791 is slightly soluble in DMSO and dimethyl formamide. It is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

PF-00489791 is an inhibitor of phosphodiesterase 5A (PDE5A; IC₅₀ = 1.9 nM).¹ It does not cross the blood-brain barrier in rats. PF-00489791 (10 mg/kg) improves motor scores in a rat model of stroke induced by middle cerebral artery occlusion (MCAO).

Reference

1. Menniti, F.S., Ren, J., Sietsma, D.K., *et al.* A non-brain penetrant PDE5A inhibitor improves functional recovery after stroke in rats. *Restor. Neurol. Neurosci.* **30(4)**, 283-289 (2012).

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, 03/15/2024

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