

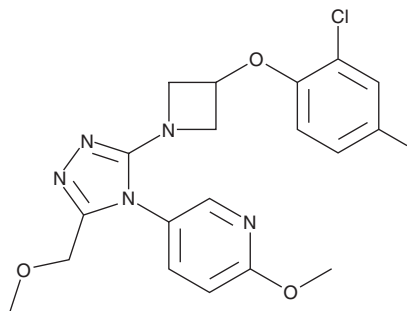
PRODUCT INFORMATION



PF-3274167

Item No. 26189

CAS Registry No.: 900510-03-4
Formal Name: 5-[3-[3-(2-chloro-4-fluorophenoxy)-1-azetidiny]-5-(methoxymethyl)-4H-1,2,4-triazol-4-yl]-2-methoxy-pyridine
MF: C₁₉H₁₉ClFN₅O₃
FW: 419.8
Purity: ≥98%
UV/Vis.: λ_{max}: 221 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-3274167 is supplied as a crystalline solid. A stock solution may be made by dissolving the PF-3274167 in the solvent of choice, which should be purged with an inert gas. PF-3274167 is soluble in the organic solvent DMSO at a concentration of approximately 18 mg/ml.

Description

PF-3274167 is an oxytocin receptor antagonist ($K_i = 9.5 \text{ nM}$).¹ It is selective for the oxytocin receptor over the vasopressin receptors V_{1A} and V_2 (K_i s = 1,120 and 10,000 nM, respectively).

Reference

1. Brown, A., Brown, T.B., Calabrese, A., *et al.* Triazole oxytocin antagonists: Identification of an aryloxyazetidiny replacement for a biaryl substituent. *Bioorg. Med. Chem. Lett.* **20(2)**, 516-520 (2010).

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, 12/19/2022

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