

PRODUCT INFORMATION



BI-639667

Item No. 41523

CAS Registry No.: 1295298-26-8

Formal Name: 1-(4-fluorophenyl)-N-[1-[2-(methylsulfonyl)-4-pyridinyl]cyclopropyl]-1H-pyrazolo[3,4-c]pyridine-4-carboxamide

MF: C₂₂H₁₈FN₅O₃S

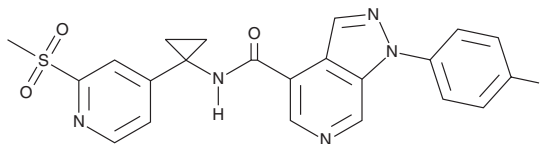
FW: 451.5

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

BI-639667 is supplied as a solid. A stock solution may be made by dissolving the BI-639667 in the solvent of choice, which should be purged with an inert gas. BI-639667 is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in acetonitrile.

Description

BI-639667 is a chemokine (C-C motif) receptor 1 (CCR1) antagonist.¹ It inhibits calcium flux induced by chemokine (C-C motif) ligand 3 (CCL3) in cells expressing CCR1 (IC₅₀ = 2 nM). BI-639667 also inhibits RANTES-induced chemotaxis in THP-1 cells (IC₅₀ = 2 nM).

Reference

1. Harcken, C., Kuzmich, D., Cook, B., *et al.* Identification of novel azaindazole CCR1 antagonist clinical candidates. *Bioorg. Med. Chem. Lett.* **29**(3), 441-448 (2019).

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, 07/30/2024

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