

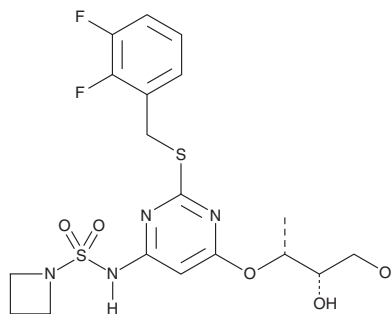
PRODUCT INFORMATION



AZD 5069

Item No. 28297

CAS Registry No.: 878385-84-3
Formal Name: N-[2-[[[(2,3-difluorophenyl)methyl]thio]-6-[[[(1R,2S)-2,3-dihydroxy-1-methylpropoxy]-4-pyrimidinyl]-1-azetidinesulfonamide
MF: C₁₈H₂₂F₂N₄O₅S₂
FW: 476.5
Purity: ≥98%
UV/Vis.: λ_{max}: 254 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

AZD 5069 is supplied as a crystalline solid. A stock solution may be made by dissolving the AZD 5069 in the solvent of choice, which should be purged with an inert gas. AZD 5069 is soluble in DMSO.

Description

AZD 5069 is an antagonist of chemokine receptor 2 (CXCR2; IC₅₀ = 0.79 nM in a radioligand binding assay).¹ It inhibits CXCL1-induced CD11b expression in and chemotaxis of isolated human neutrophils (pA₂s = 6.9 and 9.1, respectively). AZD 5069 (3 and 10 μmol/kg) inhibits bronchoalveolar lavage fluid (BALF) neutrophil influx in a rat model of LPS-induced lung neutrophilia and non-allergic airway inflammation.

Reference

1. Nicholls, D.J., Wiley, K., Dainty, I., *et al.* Pharmacological characterization of AZD5069, a slowly reversible CXCR2 chemokine receptor 2 antagonist. *J. Pharmacol. Exp. Ther.* **353**(2), 340-350 (2015).

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, 11/22/2022

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