

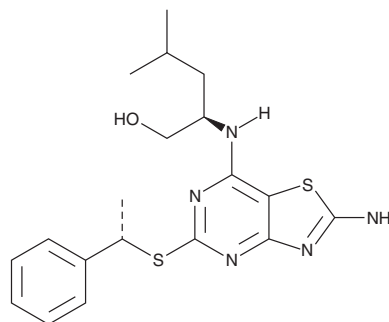
PRODUCT INFORMATION



AZD 8797

Item No. 28419

CAS Registry No.: 911715-90-7
Formal Name: (2R)-2-[[2-amino-5-[[[(1S)-1-phenylethyl]thio]thiazolo[4,5-d]pyrimidin-7-yl]amino]-4-methyl-1-pentanol
MF: C₁₉H₂₅N₅OS₂
FW: 403.6
Purity: ≥98%
UV/Vis.: λ_{max}: 246 nm
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

AZD 8797 is supplied as a solid. A stock solution may be made by dissolving the AZD 8797 in the solvent of choice, which should be purged with an inert gas. AZD 8797 is soluble in DMSO.

Description

AZD 8797 is a chemokine (C-X₃-C motif) receptor 1 (CX₃CR1) antagonist (K_i = 3.9 nM). It is selective for CX₃CR1 over CX₃CR2 (K_i = 2,800 nM).¹ AZD 8797 (0.1-10 μM) inhibits [³⁵S]GTPγS accumulation induced by chemokine (C-X₃-C) ligand 1 (CX₃CL1) in CHO cell membranes expressing human CX₃CR1.² It inhibits adhesion of RBMI-8226 cells to CX₃CL1 (IC₅₀ = 5.8 nM). AZD 8797 (39 and 78 μmol/kg per day) reduces spinal cord inflammation and demyelination, as well as the incidence of disease relapse, in a rat model of myelin oligodendrocyte glycoprotein-induced experimental autoimmune encephalomyelitis (EAE).³

References

1. Shima, H., Matsumoto, M., Ishigami, Y., *et al.* S-Adenosylmethionine synthesis is regulated by selective N⁶-adenosine methylation and mRNA degradation involving METTL16 and YTHDC1. *Cell Rep.* **21(12)**, 3354-3363 (2017).
2. Cederblad, L., Rosengren, B., Ryberg, E., *et al.* AZD8797 is an allosteric non-competitive modulator of the human CX₃CR1 receptor. *Biochem. J.* **473(5)**, 641-649 (2016).
3. Wollberg, A.R., Ericsson-Dahlstrand, A., Jureus, A., *et al.* Pharmacological inhibition of the chemokine receptor CX₃CR1 attenuates disease in a chronic-relapsing rat model for multiple sclerosis. *PNAS* **111(14)**, 5409-5414 (2014).

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, 06/28/2023

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