

PRODUCT INFORMATION



CXCR3 Antagonist 6c

Item No. 29731

CAS Registry No.: 870998-13-3

Formal Name: 4-[2-[(3-chlorobenzoyl)amino]-4-[[[2-(2,4-dichlorophenyl)ethyl]amino]carbonyl]phenyl]-N-ethylhexahydro-1H-1,4-diazepine-1-carboxamide

MF: C₃₀H₃₂Cl₃N₅O₃

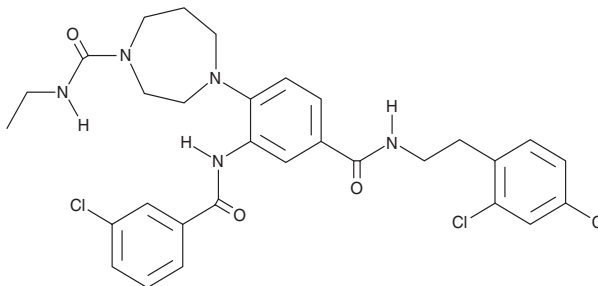
FW: 617.0

Purity: ≥98%

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

CXCR3 antagonist 6c is supplied as a crystalline solid. A stock solution may be made by dissolving the CXCR3 antagonist 6c in the solvent of choice, which should be purged with an inert gas. CXCR3 antagonist 6c is soluble in the organic solvent DMSO at a concentration of approximately 100 mM.

Description

CXCR3 antagonist 6c is an antagonist of chemokine (C-X-C motif) receptor 3 (CXCR3).¹ It inhibits calcium mobilization induced by chemokine (C-X-C motif) ligand 11 (CXCL11) in HEK293 cells expressing the human receptor (IC₅₀ = 0.06 μM). It is selective for CXCR3 over a panel of 14 human G protein-coupled receptors at 10 μM. CXCR3 antagonist 6c inhibits CXCR3-mediated migration of isolated human T cells (IC₅₀ = ~100 nM).

Reference

1. Cole, A.G., Stroke, I.L., Brescia, M.-R., *et al.* Identification and initial evaluation of 4-N-aryl-[1,4]diazepane ureas as potent CXCR3 antagonists. *Bioorg. Med. Chem. Lett.* **16(1)**, 200-203 (2006).

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, 11/30/2022

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