

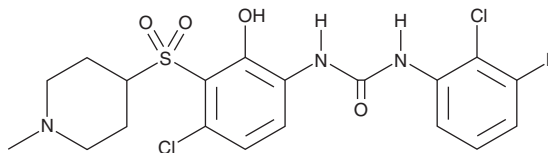
PRODUCT INFORMATION



CXCR2-IN-1

Item No. 28426

CAS Registry No.: 1873376-49-8
Formal Name: N-(2-chloro-3-fluorophenyl)-N'-[4-chloro-2-hydroxy-3-[(1-methyl-4-piperidyl)sulfonyl]phenyl]-urea
Synonym: CXCR2 Inhibitor 1
MF: C₁₉H₂₀Cl₂FN₃O₄S
FW: 476.3
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

CXCR2-IN-1 is supplied as a solid. A stock solution may be made by dissolving the CXCR2-IN-1 in the solvent of choice, which should be purged with an inert gas. CXCR2-IN-1 is soluble in DMSO.

Description

CXCR2-IN-1 is a chemokine (C-X-C motif) receptor 2 (CXCR2) antagonist (IC₅₀ = 0.501 nM in a reporter assay).¹ It inhibits chemotaxis of isolated human neutrophils induced by chemokine (C-X-C motif) ligand 1 (CXCL1; IC₅₀ = 79.4 nM). *In vivo*, CXCR2-IN-1 (100 mg/kg) inhibits forebrain lesion formation in a mouse model of cuprizone-induced demyelination.

Reference

- Xu, H., Lu, H., Xu, Z., *et al.* Discovery of CNS penetrant CXCR2 antagonists for the potential treatment of CNS demyelinating disorders. *Med. Chem. Lett.* **7**(4), 397-402 (2016).

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, 03/25/2024

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