

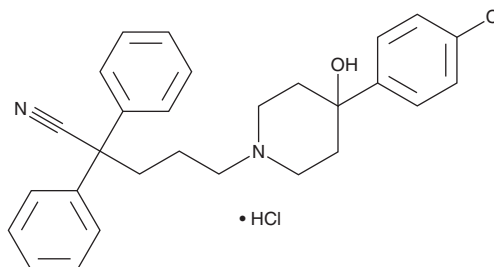
PRODUCT INFORMATION



BX-513 (hydrochloride)

Item No. 42733

CAS Registry No.: 1216540-18-9
Formal Name: 4-(4-chlorophenyl)-4-hydroxy- α,α -diphenyl-1-piperidinepentanenitrile, monohydrochloride
Synonym: VUF 2274
MF: $C_{28}H_{29}ClN_2O \cdot HCl$
FW: 481.5
Purity: $\geq 98\%$
Supplied as: A solid
Storage: $-20^\circ C$
Stability: ≥ 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

BX-513 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the BX-513 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. BX-513 (hydrochloride) is slightly soluble (0.1-1 mg/ml) in ethanol and sparingly soluble (1-10 mg/ml) in DMSO.

Description

BX-513 is an antagonist of chemokine (C-C motif) receptor 1 (CCR1; $K_i = 0.04 \mu M$).¹ It is selective for CCR1 over 15 other receptors (K_i s = $2 \rightarrow 10 \mu M$). BX-513 inhibits increases in acidification induced by chemokine (C-C motif) ligand 18 (CCL18) in THP-1 monocytes in a concentration-dependent manner. It inhibits RANTES- and CCL18-induced chemotaxis in isolated peripheral blood mononuclear cells (PBMCs) when used at concentrations of 10 or 100 nM.

Reference

1. Hesselgesser, J., Ng, H.P., Liang, M., *et al.* Identification and characterization of small molecule functional antagonists of the CCR1 chemokine receptor. *J. Biol. Chem.* **273**(25), 15687-15692 (1998).

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, 02/11/2025

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