

PRODUCT INFORMATION



CCR2 Antagonist

Item No. 19542

CAS Registry No.: 445479-97-0

Formal Name: *rel*-2-[[[(1-methylethyl)amino]carbonyl]amino]-N-[2-[[[(1R,2S)-2-[[4-(methylthio)benzoyl]amino]cyclohexyl]amino]-2-oxoethyl]-5-(trifluoromethyl)-benzamide

Synonym: BMS CCR2-22

MF: $C_{28}H_{34}F_3N_5O_4S$

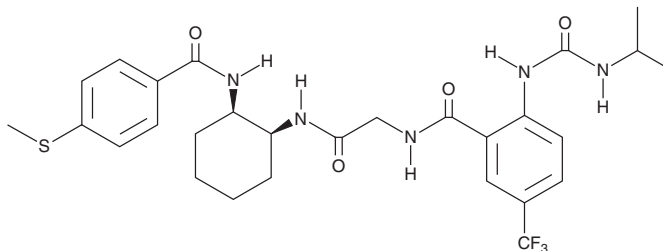
FW: 593.7

Purity: $\geq 98\%$

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

CCR2 antagonist is supplied as a solid. A stock solution may be made by dissolving the CCR2 antagonist in the solvent of choice, which should be purged with an inert gas. CCR2 antagonist is soluble (≥ 10 mg/ml) in DMSO.

Description

CCR2 antagonist is an antagonist of chemokine (C-C motif) receptor 2 (CCR2).¹ It binds to CCR2 in a radioligand binding assay using peripheral blood mononuclear cells (PBMCs; $IC_{50} = 5.1$ nM). CCR2 antagonist inhibits chemotaxis induced by chemokine (C-C motif) ligand 2 (CCL2) in PBMCs ($IC_{50} = 1$ nM). *In vivo*, CCR2 antagonist (2 mg/kg) reduces hepatic Ly6C⁺ monocyte recruitment and viral clearance in mice infected with hepatitis B virus (HBV).²

References

- Cherney, R.J., Mo, R., Meyer, D.T., *et al.* Discovery of disubstituted cyclohexanes as a new class of CC chemokine receptor 2 antagonists. *J. Med. Chem.* **51**(4), 721-724 (2008).
- Wu, L.L., Peng, W.H., Wu, H.L., *et al.* Lymphocyte antigen 6 complex, locus C⁺ monocytes and Kupffer cells orchestrate liver immune responses against hepatitis B virus in mice. *Hepatology* **69**(6), 2364-2380 (2019).

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, 10/07/2024

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