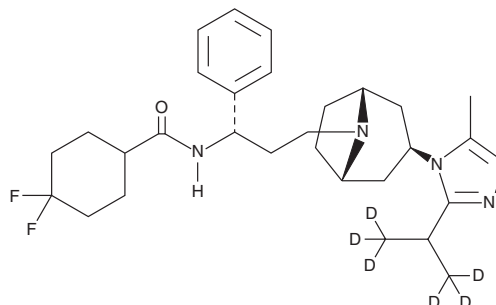


PRODUCT INFORMATION



Maraviroc-d₆ Item No. 28767

CAS Registry No.: 1033699-22-7
Formal Name: 4,4-difluoro-N-[(1S)-3-[(3-exo)-3-[3-methyl-5-(1-(methyl-d₃)ethyl-2,2,2-d₃)-4H-1,2,4-triazol-4-yl]-8-azabicyclo[3.2.1]oct-8-yl]-1-phenylpropyl]-cyclohexanecarboxamide
MF: C₂₉H₃₅D₆F₂N₅O
FW: 519.7
Chemical Purity: ≥98% (Maraviroc)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₆); ≤1% d₀
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

Maraviroc-d₆ is intended for use as an internal standard for the quantification of maraviroc (Item No. 14641) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Maraviroc-d₆ is supplied as a solid. A stock solution may be made by dissolving the maraviroc-d₆ in the solvent of choice, which should be purged with an inert gas. Maraviroc-d₆ is slightly soluble in chloroform.

Description

Maraviroc is an antagonist of the chemokine (C-C motif) receptor 5 (CCR5; K_i = 0.24 nM for the rhesus monkey recombinant receptor).¹ It inhibits binding of chemokine (C-C motif) ligand 3 (CCL3), CCL4, and CCL5 to HEK293 cell membranes expressing CCR5 (IC₅₀s = 3.3, 7.2, and 5.2 nM, respectively) and inhibits HIV-1 binding to CCR5 via glycoprotein 120 (gp120) and gp160 (IC₅₀s = 11 and 0.22 nM, respectively).² Maraviroc is selective for CCR5 over CCR1, -2, -3, -4, -7, and -8, and chemokine (C-X-C motif) receptor 1 (CXCR1) and CXCR2 in a panel of immunological assays for ligand-induced cell chemotaxis and ligand-receptor binding (IC₅₀s = >10 μM for all). It has antiviral activity against laboratory and clinical isolates of CCR5-tropic (IC₅₀s = 0.1-1.1 nM), but not CXCR4-tropic or dual tropic (IC₅₀s = >10 μM for all), HIV-1 in isolated human peripheral blood mononuclear cells (PBMCs). Maraviroc prevents infection upon exposure to HIV-1 in humanized RAG-hum mice when administered at a dose of 62 mg/kg.³

References

1. Yuan, Y., Arnatt, C.K., El-Hage, N., *et al.* A bivalent ligand targeting the putative mu opioid receptor and chemokine receptor CCR5 heterodimers: Binding affinity versus functional activities. *Medchemcomm.* **4**(5), 847-851 (2013).
2. Dorr, P., Westby, M., Dobbs, S., *et al.* Maraviroc (UK-427,857), a potent, orally bioavailable, and selective small-molecule inhibitor of chemokine receptor CCR5 with broad-spectrum anti-human immunodeficiency virus type 1 activity. *Antimicrob. Agents Chemother.* **49**(11), 4721-4732 (2005).
3. Neff, C.P., Ndolo, T., Tandon, A., *et al.* Oral pre-exposure prophylaxis by anti-retrovirals raltegravir and maraviroc protects against HIV-1 vaginal transmission in a humanized mouse model *PLoS One* **5**(12), e15257 (2010).

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, 03/07/2023

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