

PRODUCT INFORMATION

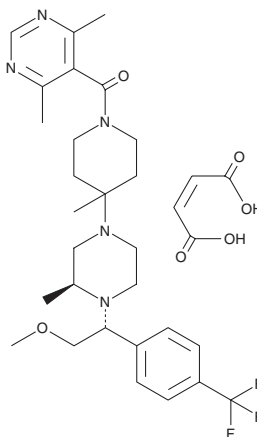


Viciviroc (maleate)

Item No. 28453

CAS Registry No.: 599179-03-0
Formal Name: 1-[(4,6-dimethyl-5-pyrimidinyl)carbonyl]-4-[(3S)-4-[(1R)-2-methoxy-1-[4-(trifluoromethyl)phenyl]ethyl]-3-methyl-1-piperazinyl]-4-methyl-piperidine, 2Z-butenedioate

Synonym: SCH 417690
MF: C₂₈H₃₈F₃N₅O₂ • C₄H₄O₄
FW: 649.7
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

Viciviroc (maleate) is supplied as a crystalline solid. A stock solution may be made by dissolving the viciviroc (maleate) in the solvent of choice, which should be purged with an inert gas. Viciviroc (maleate) is slightly soluble in chloroform, methanol, and DMSO (warmed).

Description

Viciviroc is a chemokine (C-C motif) receptor 5 (CCR5) antagonist ($K_i = 2.5$ nM).¹ It selectively inhibits CCR5 over M₁ and M₂ muscarinic acetylcholine receptors ($K_s = >10,000$ nM for both) and human-ether-a-go-go (hERG) potassium channels ($IC_{50} = 5.8$ μM).¹ Viciviroc decreases chemotaxis induced by chemokine (C-C motif) ligand 3 (CCL3) in Ba/F3 cells expressing recombinant human CCR5 with an IC_{50} value of 0.91 nM.² It inhibits the replication of 30 R5-tropic HIV-1 clinical isolates in peripheral blood mononuclear cells (PBMCs; $EC_{50s} = 0.04$ -1.4 nM) and reduces viral entry of seven drug-resistant strains of HIV-1 in U87 cells expressing CD4 and CCR5 ($EC_{50s} = 8.7$ -32.9 nM).

References

1. Tagat, J.R., McCombie, S.W., Nazareno, D., *et al.* Piperazine-based CCR5 antagonists as HIV-1 inhibitors. IV. Discovery of 1-[(4,6-dimethyl-5-pyrimidinyl)carbonyl]-4-[4-[2-methoxy-1(R)-4-(trifluoromethyl)phenyl]ethyl]-3(S)-methyl-1-piperazinyl]-4-methylpiperidine (Sch-417690/Sch-D), a potent, highly selective, and orally bioavailable CCR5 antagonist. *J. Med. Chem.* **47(10)**, 2405-2408 (2004).
2. Strizki, J.M., Tremblay, C., Xu, S., *et al.* Discovery and characterization of viciviroc (SCH 417690), a CCR5 antagonist with potent activity against human immunodeficiency virus type 1. *Antimicrob. Agents Chemother.* **49(12)**, 4911-4919 (2005).

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, 11/29/2022

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