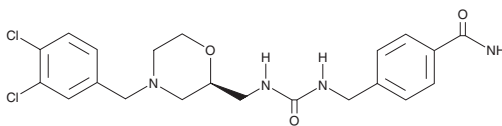


PRODUCT INFORMATION



GW 766994
Item No. 28431

CAS Registry No.: 408303-43-5
Formal Name: 4-[[[[(2S)-4-[(3,4-dichlorophenyl)methyl]-2-morpholinyl)methyl]amino]carbonyl]amino]methyl]-benzamide
Synonym: GW 994
MF: C₂₁H₂₄Cl₂N₄O₃
FW: 451.4
Purity: ≥98%
UV/Vis.: λ_{max}: 225 nm
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

GW 766994 is supplied as a solid. A stock solution may be made by dissolving the GW 766994 in the solvent of choice, which should be purged with an inert gas. GW 766994 is soluble in DMSO.

Description

GW 766994 is an antagonist of chemokine (C-C motif) receptor 3 (CCR3) with a K_i value of 13.8 nM in an assay of eosinophil chemotaxis induced by chemokine (C-C motif) ligand 11 (CCL11).¹ It decreases CCL11-induced cyclin-dependent kinase 5 (Cdk5) and tau phosphorylation and production of amyloid β (1-42) (Aβ42) in isolated mouse hippocampal neurons when used at a concentration of 10 μM.²

References

1. Neighbour, H., Boulet, L.-P., Lemiere, C., *et al.* Safety and efficacy of an oral CCR3 antagonist in patients with asthma and eosinophilic bronchitis: A randomized, placebo-controlled clinical trial. *Clin. Exp. Allergy* **44**(4), 508-516 (2014).
2. Zhu, C., Xu, B., Sun, X., *et al.* Targeting CCR3 to reduce amyloid-β production, tau hyperphosphorylation, and synaptic loss in a mouse model of Alzheimer's disease. *Mol. Neurobiol.* **54**(10), 7964-7978 (2017).

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, 09/29/2022

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