

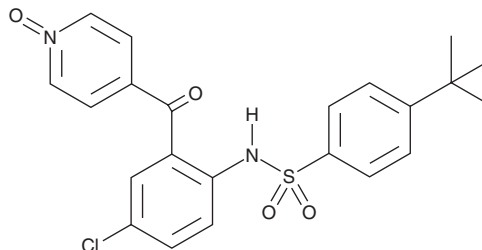
PRODUCT INFORMATION



Vercirnon

Item No. 23131

CAS Registry No.: 698394-73-9
Formal Name: N-[4-chloro-2-[(1-oxido-4-pyridinyl)carbonyl]phenyl]-4-(1,1-dimethylethyl)-benzenesulfonamide
Synonyms: CCX282-B, GSK1605786
MF: C₂₂H₂₁ClN₂O₄S
FW: 444.9
Purity: ≥98%
UV/Vis.: λ_{max}: 226, 313 nm
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

Vercirnon is supplied as a crystalline solid. A stock solution may be made by dissolving the vercirnon in the solvent of choice, which should be purged with an inert gas. Vercirnon is soluble in the organic solvent DMSO at a concentration of approximately 25 mg/ml.

Description

Vercirnon is an orally bioavailable antagonist of chemokine receptor 9 (CCR9) with an IC₅₀ value of 5.4 nM for inhibition of CCL25-induced calcium mobilization in Molt-4 cells.¹ It is selective for CCR9 over CCR1-12 and CX3CR1-7 (IC₅₀s = >10 μM for all). It also inhibits CCL25-induced Molt-4 chemotaxis in 0.1% BSA/HBSS and in 100% human AB serum (IC₅₀s = 3.5 and 33.4 nM, respectively), as well as chemotaxis of human T cells stimulated by retinoic acid (RA) and mouse and rat thymocytes. Vercirnon binds allosterically to the intracellular side of CCR9 and prevents G protein coupling.²

References

1. Walters, M.J., Wang, Y., Lai, N., *et al.* Characterization of CCX282-B, an orally bioavailable antagonist of the CCR9 chemokine receptor, for treatment of inflammatory bowel disease. *J. Pharmacol. Exp. Ther.* **335**(1), 61-69 (2010).
2. Oswald, C., Rappas, M., Kean, J., *et al.* Intracellular allosteric antagonism of the CCR9 receptor. *Nature* **540**(7633), 462-465 (2016).

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, 10/06/2022

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