

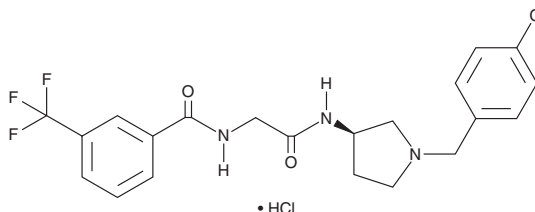
PRODUCT INFORMATION



Teijin Compound 1 (hydrochloride)

Item No. 30977

CAS Registry No.: 1313730-14-1
Formal Name: N-[2-[[[(3R)-1-[(4-chlorophenyl)methyl]-3-pyrrolidinyl]amino]-2-oxoethyl]-3-(trifluoromethyl)benzamide, monohydrochloride
MF: C₂₁H₂₁ClF₃N₃O₂ • HCl
FW: 476.3
Purity: ≥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

Teijin compound 1 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the teijin compound 1 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Teijin compound 1 (hydrochloride) is soluble in the organic solvent DMSO. It is also soluble in water. The solubility of teijin compound 1 (hydrochloride) in DMSO and water is approximately 100 and 10 mM, respectively. We do not recommend storing the aqueous solution for more than one day.

Description

Teijin compound 1 is an antagonist of chemokine (C-C motif) receptor 2b (CCR2b; IC₅₀ = 180 nM in a radioligand binding assay).^{1,2} It inhibits chemotaxis induced by chemokine (C-C motif) ligand 2 (CCL2) in THP-1 cells with an EC₅₀ value of 24 nM.¹

References

1. Moree, W.J., Kataoka, K.-i., Ramirez-Weinhouse, M.M., *et al.* Small molecule antagonists of the CCR2b receptor. Part 2: Discovery process and initial structure-activity relationships of diamine derivatives. *Bioorg. Med. Chem. Lett.* **14**(21), 5413-5416 (2004).
2. Hall, S.E., Mao, A., Nicolaidou, V., *et al.* Elucidation of binding sites of dual antagonists in the human chemokine receptors CCR2 and CCR5. *Mol. Pharmacol.* **75**(6), 1325-1336 (2009).

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, 11/04/2022

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