

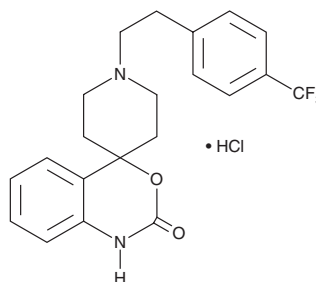
PRODUCT INFORMATION



RS 102895 (hydrochloride)

Item No. 18207

CAS Registry No.: 1173022-16-6
Formal Name: 1'-[2-[4-(trifluoromethyl)phenyl]ethyl]-spiro[4H-3,1-benzoxazine-4,4'-piperidin]-2(1H)-one, monohydrochloride
MF: C₂₁H₂₁F₃N₂O₂ • HCl
FW: 426.9
Purity: ≥98%
UV/Vis.: λ_{max}: 204, 279 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

RS 102895 (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the RS 102895 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. RS 102895 (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of RS 102895 (hydrochloride) in these solvents is approximately 2, 30, and 25 mg/ml, respectively.

RS 102895 (hydrochloride) is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, RS 102895 (hydrochloride) should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. RS 102895 (hydrochloride) has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

The chemokine CCL2, also known as monocyte chemoattractant protein-1 (MCP-1), stimulates leukocyte chemotaxis to sites of inflammation via signaling through the MCP-1 receptor, CCR2. RS 102895 is a spiro-piperidine compound that acts as a CCR2 antagonist (IC₅₀ = 0.36 μM).¹ It inhibits the related CCR1 receptor with an IC₅₀ value of 17.8 μM and inhibits adrenergic α_{1a}, α_{1d}, and 5HT_{1A} receptors with IC₅₀ values of 0.13, 0.32, and 47 μM, respectively.¹ RS 102895 prevents chemotaxis of THP-1 cells (IC₅₀ = 1.7 μM) when MCP-1 is presented as a chemoattractant.¹

Reference

1. Mirzadegan, T., Diehl, F., Ebi, B., *et al.* Identification of the binding site for a novel class of CCR2b chemokine receptor antagonists: Binding to a common chemokine receptor motif within the helical bundle. *J. Biol. Chem.* **275**(33), 25562-25571 (2000).

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, 11/01/2022

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