

PRODUCT INFORMATION

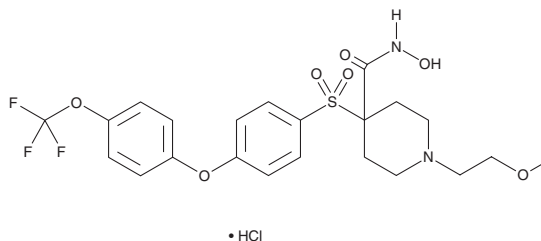


SD 2590 (hydrochloride)

Item No. 27680

CAS Registry No.: 226395-93-3
Formal Name: N-hydroxy-1-(2-methoxyethyl)-4-[[4-[4-(trifluoromethoxy)phenoxy]phenyl]sulfonyl]-4-piperidinecarboxamide, monohydrochloride

Synonym: SC-78080
MF: C₂₂H₂₅F₃N₂O₇S • HCl
FW: 555.0
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

SD 2590 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the SD 2590 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. SD 2590 (hydrochloride) is soluble in organic solvents such as DMSO. It is also soluble in water. The solubility of SD 2590 (hydrochloride) in DMSO and water is approximately 100 mM. We do not recommend storing the aqueous solution for more than one day.

Description

SD 2590 is an inhibitor of matrix metalloproteinase-2 (MMP-2), -3, -8, -9, and -13 (IC₅₀s = <0.1, 28.7, 1.7, 0.18, and <0.1 nM, respectively).¹ It is selective for these MMPs over MMP-1 and MMP-7 (IC₅₀s = >10,000 and 7,000 nM, respectively). SD 2590 (10 mg/kg per day) inhibits dilation of the left ventricle in rats following myocardial infarction (MI) compared to MI vehicle control animals.

Reference

1. Becker, D.P., Barta, T.E., Bedell, L.J., *et al.* Orally active MMP-1 sparing α -tetrahydropyranyl and α -piperidinyl sulfone matrix metalloproteinase (MMP) inhibitors with efficacy in cancer, arthritis, and cardiovascular disease. *J. Med. Chem.* **53**(18), 6653-6680 (2010).

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, 12/01/2022

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