

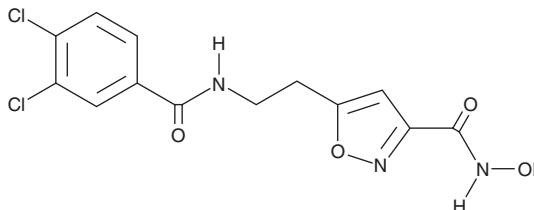
PRODUCT INFORMATION



SS-208

Item No. 29128

CAS Registry No.: 2245942-72-5
Formal Name: 5-[2-[(3,4-dichlorobenzoyl)amino]ethyl]-N-hydroxy-3-isoxazolecarboxamide
MF: C₁₃H₁₁Cl₂N₃O₄
FW: 344.2
Purity: ≥98%
UV/Vis.: λ_{max}: 237 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

SS-208 is supplied as a crystalline solid. A stock solution may be made by dissolving the SS-208 in the solvent of choice, which should be purged with an inert gas. SS-208 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of SS-208 in ethanol is approximately 30 mg/ml and approximately 1 mg/ml in DMSO and DMF.

SS-208 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, SS-208 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. SS-208 has a solubility of approximately 0.3 mg/ml in a 1:2 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

SS-208 is an inhibitor of histone deacetylase 6 (HDAC6; IC₅₀ = 12 nM).¹ It is selective for HDAC6 over HDAC1, -4, -5, -7, -8, -9, and -11 (IC₅₀s = 1.39, 19.5, 6.91, 8.34, 1.23, 38.2, and 5.12 μM, respectively). SS-208 (0.1-25 μM) inhibits HDAC6 in, but does not induce cell death of, human PC3 prostate, 5637 and T24 bladder, and murine SM1 melanoma cells *in vitro*. It decreases protein levels of programmed death ligand 1 (PD-L1) in SM1 cells when used at a concentration of 5 μM. SS-208 (25 mg/kg, i.p.) reduces tumor growth and increases the number of intratumoral CD8⁺, CD4⁺, and natural killer (NK) T cells in an SM1 murine melanoma model.

Reference

1. Shen, S., Hadley, M., Ustinova, K., *et al.* Discovery of a new isoxazole-3-hydroxamate-based histone deacetylase 6 inhibitor SS-208 with antitumor activity in syngeneic melanoma mouse models. *J. Med. Chem.* **62**(18), 8557-8577 (2019).

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, 03/05/2020

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