

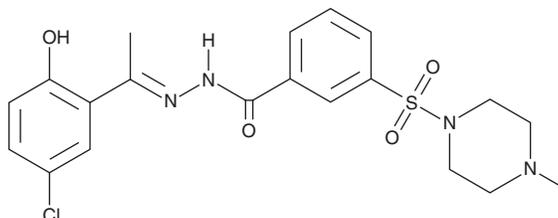
PRODUCT INFORMATION



SP-2577

Item No. 43468

CAS Registry No.: 1423715-37-0
Formal Name: 3-[(4-methyl-1-piperazinyl)sulfonyl]-benzoic acid, (2E)-2-[1-(5-chloro-2-hydroxyphenyl)ethylidene]hydrazide
Synonym: Seclidemstat
MF: C₂₀H₂₃ClN₄O₄S
FW: 450.9
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

SP-2577 is supplied as a solid. A stock solution may be made by dissolving the SP-2577 in the solvent of choice, which should be purged with an inert gas. SP-2577 is soluble (≥10 mg/ml) in DMSO. SP-2577 is slightly soluble (0.1-1 mg/ml) in ethanol.

Description

SP-2577 is an inhibitor of lysine-specific demethylase 1 (LSD1; IC₅₀ = 13 nM).¹ It induces cell death in a panel of 12 ovarian cancer cell lines expressing mutant switch/sucrose nonfermentable (SWI/SNF; IC₅₀s = 0.013-2.819 μM). SP-2577 increases secretion of IFN-γ, IFN-β, IL-1β, IL-2, and IL-6 in COV434 ovarian cancer cells. It decreases the viability of primary natural killer T (NKT) cells and T cells and reduces the mitochondrial oxygen consumption rate (OCR) in NKT cells when used at a concentration of 5 μM.² SP-2577 (5 μM) induces differentiation of primary rat neural stem cells into astrocytes.³ It increases infiltration of human peripheral blood mononuclear cells (PBMCs) into isolated COV434 tumor organoids in a migration and invasion assay when used at a concentration of 3 μM.¹

References

1. Soldi, R., Halder, T.G., Weston, A., *et al.* The novel reversible LSD1 inhibitor SP-2577 promotes anti-tumor immunity in SWI/SNF complex mutated ovarian cancer. *PLoS One* **15**(7), e0235705 (2020).
2. Bailey, C.P., Figueroa, M., Gangadharan, A., *et al.* Scaffolding LSD1 inhibitors impair NK cell metabolism and cytotoxic function through depletion of glutathione. *Front. Immunol.* **11**, 2196 (2020).
3. Kim, S. and Kim, H.-J. Histone lysine demethylase 1A inhibitors, seclidemstat and tranilcypropridine, induce astrocytogenesis in rat neural stem cells. *Biochem. Biophys. Res. Commun.* **750**, 151330 (2025).

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, 06/17/2025

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