

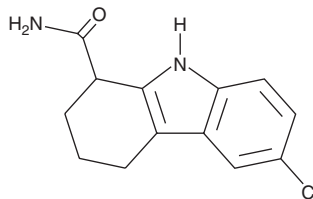
PRODUCT INFORMATION



(±)-EX-527

Item No. 10009798

CAS Registry No.: 49843-98-3
Formal Name: 6-chloro-2,3,4,9-tetrahydro-1H-carbazole-1-carboxamide
MF: C₁₃H₁₃ClN₂O
FW: 248.7
Purity: ≥98%
UV/Vis.: λ_{max}: 233, 291 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

(±)-EX-527 is supplied as a crystalline solid. A stock solution may be made by dissolving the (±)-EX-527 in the solvent of choice, which should be purged with an inert gas. (±)-EX-527 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of (±)-EX-527 in ethanol is approximately 5 mg/ml and approximately 20 mg/ml in DMSO and DMF.

(±)-EX-527 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, (±)-EX-527 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. (±)-EX-527 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

EX-527 is an inhibitor of sirtuin 1 (SIRT1; IC₅₀ = 0.098 μM).¹ It is selective for SIRT1 over SIRT2, and SIRT3 (IC₅₀s = 19.6 and 48.7 μM, respectively) and the cytochrome P450 (CYP) isoforms CYP3A4, CYP2D6, CYP2C9, CYP2C19, and CYP1A2 at 1 μM. EX-527 (50 μM) induces cell cycle arrest at the G₁ phase in MCF-7 cells.²

References

1. Napper, A.D., Hixon, J., McDonagh, T., *et al.* Discovery of indoles as potent and selective inhibitors of the deacetylase SIRT1. *J. Med. Chem.* **48(25)**, 8045-8054 (2005).
2. Peck, B., Chen, C.-Y., Ho, K.-K., *et al.* SIRT inhibitors induce cell death and p53 acetylation through targeting both SIRT1 and SIRT2. *Mol. Cancer Ther.* **9(4)**, 844-855 (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, 10/10/2022

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