

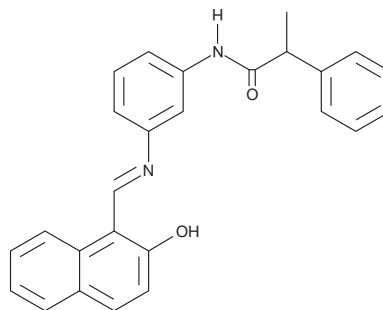
PRODUCT INFORMATION



Salermide

Item No. 13178

CAS Registry No.: 1105698-15-4
Formal Name: N-[3-[[[2-hydroxy-1-naphthalenyl)methylene]amino]phenyl]-a-methylbenzeneacetamide
MF: C₂₆H₂₂N₂O₂
FW: 394.5
Purity: ≥98%
UV/Vis.: λ_{max}: 207, 233, 318, 365, 440, 459 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

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

Salermide is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, salermide should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Salermide has a solubility of approximately 0.25 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

Salermide is an inhibitor of SIRT1 and SIRT2, causing tumor-specific apoptotic cell death.¹ In MOLT4 leukemia cells, salermide causes 90% apoptosis within 72 hours (IC₅₀ ~20 μM) by reactivating proapoptotic genes that are repressed by SIRT1.¹

Reference

1. Yamamoto, H., Schoonjans, K., and Auwerx, J. Sirtuin functions in health and disease. *Mol. Endocrinol.* **21(8)**, 1745-1755 (2007).
2. Milne, J.C. and Denu, J.M. The sirtuin family: Therapeutic targets to treat diseases of aging. *Curr. Opin. Chem. Biol.* **12(1)**, 11-17 (2008).
3. Lara, E., Mai, A., Calvanese, V., et al. Salermide, a sirtuin inhibitor with a strong cancer-specific proapoptotic effect. *Oncogene* **28(6)**, 781-791 (2008).

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, 03/11/2024

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