

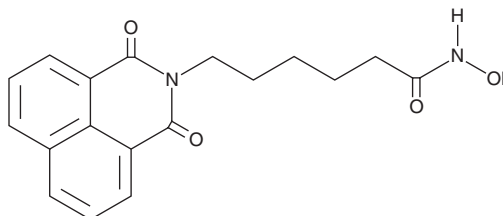
PRODUCT INFORMATION



Scriptaid

Item No. 10572

CAS Registry No.: 287383-59-9
Formal Name: N-hydroxy-1,3-dioxo-1H-benz[de]isoquinoline-2(3H)-hexanamide
Synonym: GCK 1026
MF: C₁₈H₁₈N₂O₄
FW: 326.4
Purity: ≥98%
UV/Vis.: λ_{max}: 234, 333 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

Scriptaid is supplied as a crystalline solid. A stock solution may be made by dissolving the scriptaid in the solvent of choice, which should be purged with an inert gas. Scriptaid is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of scriptaid in these solvents is approximately 2.5 and 3 mg/ml, respectively.

Scriptaid is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, scriptaid should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Scriptaid has a solubility of approximately 0.1 mg/ml in a 1:9 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Scriptaid is a histone deacetylase inhibitor that has an optimal concentration of 6-8 μM in a cell-based assay, is less toxic than trichostatin A, and works in a wide variety of biological systems.¹ It induces cell cycle arrest in colon cancer cells in culture and inhibits tumor growth *in vitro* and *in vivo*.^{2,3} Scriptaid also facilitates the cloning of inbred mouse strains produced by somatic cell nuclear transfer.⁴

References

1. Su, G.H., Sohn, T.A., Ryu, B., *et al.* A novel histone deacetylase inhibitor identified by high-throughput transcriptional screening of a compound library. *Cancer Res.* **60(12)**, 3137-3142 (2000).
2. Keen, J.C., Yan, L., Mack, K.M., *et al.* A novel histone deacetylase inhibitor, scriptaid, enhances expression of functional estrogen receptor α (ER) in ER negative human breast cancer cells in combination with 5-aza 2'-deoxycytidine. *Breast Cancer Res. Treat.* **81(3)**, 177-186 (2003).
3. Lee, E.J., Lee, B.B., Kim, S.J., *et al.* Histone deacetylase inhibitor scriptaid induces cell cycle arrest and epigenetic change in colon cancer cells. *Int. J. Oncol.* **33(4)**, 767-776 (2008).
4. Van Thuan, N., Bui, H.T., Kim, J.H., *et al.* The histone deacetylase inhibitor scriptaid enhances nascent mRNA production and rescues full-term development in cloned inbred mice. *Reproduction* **138(2)**, 309-317 (2009).

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, 11/17/2022

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