

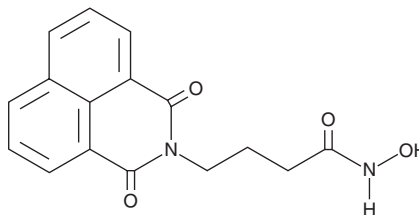
PRODUCT INFORMATION



Nullscript

Item No. 16433

CAS Registry No.: 300816-11-9
Formal Name: N-hydroxy-1,3-dioxo-1H-benz[de]isoquinoline-2(3H)-butanamide
MF: C₁₆H₁₄N₂O₄
FW: 298.3
Purity: ≥98%
UV/Vis.: λ_{max}: 212, 234, 332, 344 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

Nullscript is supplied as a crystalline solid. A stock solution may be made by dissolving the nullscript in the solvent of choice, which should be purged with an inert gas. Nullscript is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of nullscript in these solvents is approximately 2 mg/ml.

Nullscript is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, nullscript should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Nullscript 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 (Item No. 10572) is a histone deacetylase (HDAC) 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 Nullscript is an analog of scriptaid that lacks the HDAC inhibitory effects of scriptaid.¹ It is used as a negative control in experiments involving scriptaid.¹

Reference

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).

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, 01/03/2023

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