

Scriptaid (GCK 1026)

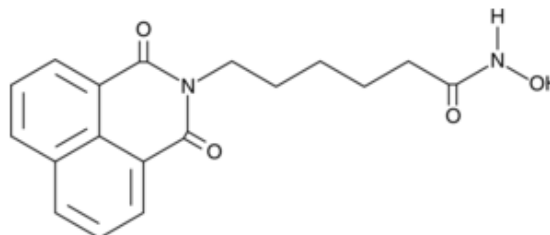
Catalog #: 27200

Lot #: 120203

Size: 1 mg

Structure:

CAS Registry #: 287383-59-9



Purity: ≥98%

Chemical Formula: C₁₈H₁₈N₂O₄

Molecular Weight: 326.4

Description: Scriptaid, also known as GCK 1026, is a histone deacetylase inhibitor with optimal concentration of 6-8 μM in a cell based assay. Scriptaid exhibits lower toxicity than trichostatin A. Scriptaid inhibits tumor growth *in vitro* and *in vivo*. It also induces cell cycle arrest in colon cancer cells in culture.

Appearance: White to off-white crystalline solid

Solubility: Solubility in DMSO and DMF is 2.5 mg/ml. Sparingly soluble in aqueous buffers. For maximum solubility, dissolve first in DMF and then dilute with aqueous buffer of choice. Do not store in aqueous solution for more than one day.

Biological Activity: Inhibits HDAC1 and HDAC3 with an IC₅₀ value ~0.6 μM. Also inhibits HDAC8 with an IC₅₀ ~ 1 μM.

Storage/Stability: Store at or below -20°C.

Quality Control: The purity was determined by HPLC analysis.

References:

1. Su, G.H., *et al.*, *Cancer Res.* 2010; **60**: 3137-3142.
2. Keen, J.C., *et al.*, *Cancer Res. Treat.* 2003; **81**: 177-186.
3. Lee, E.J., *et al.*, *Int. J. Oncol.* 2008; **33**: 767-776.