

Nullscript

HDAC inhibitor

Negative control compound for the histone deacetylase (HDAC) inhibitor scriptaid (Prod. No. BML-GR326).

Citations: 2

[View Online »](#)

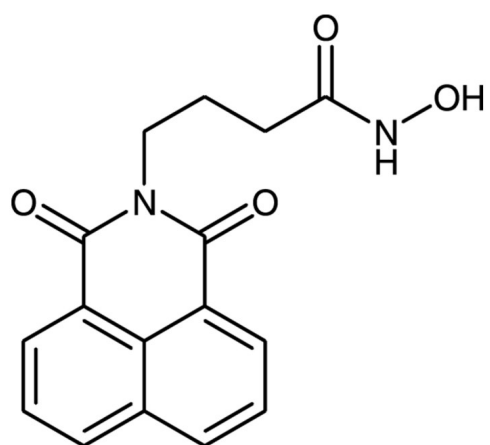
Ordering Information

[Order Online »](#)

BML-GR327-0001	1mg
----------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored, as supplied, at -20°C. Stock solutions are stable for up to 3 months at -20°C.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	4-(1,3-Dioxo-1H,3H-benzo[de]isoquinolin-2-yl)-N-hydroxybutanamide
Appearance	Yellow to off-white solid.
CAS	300816-11-9
Couple Target	HDAC
Couple Type	Inhibitor
Formula	$C_{16}H_{14}N_2O_4$
MW	298.3
Purity	≥97% (TLC)
Solubility	Soluble in DMSO (4mg/ml) or ethanol (0.3mg/ml).
Source	Synthetic.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com