

Scriptaid

HDAC inhibitor

Inhibitor of histone deacetylase (HDAC) (HDAC1 and 3: IC₅₀~0.6μM; HDAC8: IC₅₀~1μM) with lower toxicity than trichostatin A (Prod. No. BML-GR309). Inhibits tumor growth *in vitro* and *in vivo* and, in conjunction with 5-aza-2'-deoxycytidine (Prod. No. ALX-480-096), enhances expression of estrogen receptor α. Optimal concentration 2-2.5μg/ml (6-8μM). A negative control is available (Prod. No. BML-GR327).

Citations: 8

[View Online »](#)

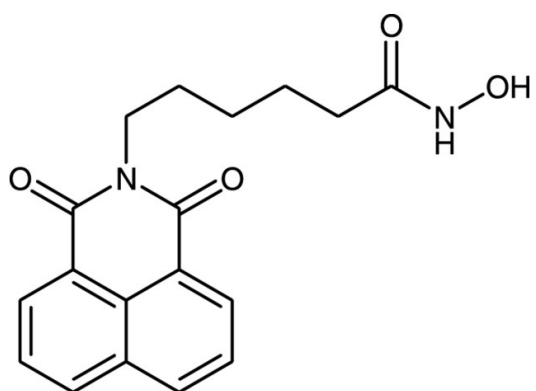
Ordering Information

[Order Online »](#)

BML-GR326-0001	1mg
BML-GR326-0005	5mg

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 air. Store under inert gas
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	6-(1,3-Dioxo-1H,3H-benzo[de]isoquinolin-2-yl)-N-hydroxyhexanamide
Appearance	Yellow solid.
CAS	287383-59-9
Couple Target	HDAC, HDAC1, HDAC3, HDAC8
Couple Type	Inhibitor
Formula	$C_{18}H_{18}N_2O_4$
Identity	Determined by NMR.
MW	326.4
Purity	≥97% (HPLC)
Solubility	Soluble in DMSO (4mg/ml), ethanol (0.3mg/ml) or methanol.

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-fr@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