

Niltubacin

HDAC inhibitor

Negative control for Tubacin (tubulin acetylation inducer), a highly potent and selective, reversible, cell-permeable inhibitor of HDAC6. Concentration of niltubacin in cell culture experiments typically ranges from 2-20µM.

Citations: 5

[View Online »](#)

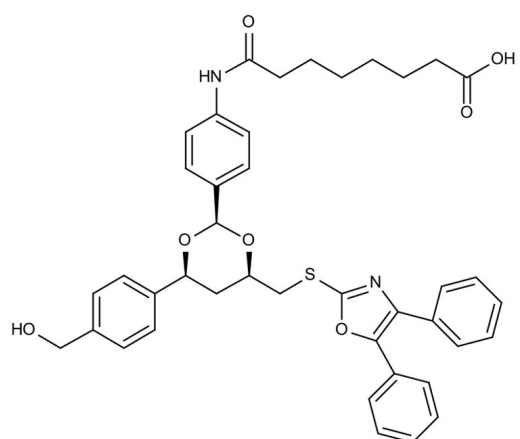
Ordering Information

[Order Online »](#)

BML-GR363-0100	100µg
BML-GR363-0500	500µg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least one year after receipt when stored at -80°C.
Short Term Storage	-20°C
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	7-(4-((2R,4R,6S)-4-((4,5-dipenyloxazol-2-ylthio)methyl)-6-(4-(hydroxymethyl)phenyl)-1,3-dioxan-2-yl)phenylcarbamoyl)heptanoic acid
Appearance	White solid.
Couple Target	HDAC, HDAC6
Couple Type	Inhibitor
Formula	C ₄₁ H ₄₂ N ₂ O ₇ S
MW	706.9
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (6 mg/ml).
Source	Synthetic.



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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