

Tubacin

HDAC inhibitor

Tubacin (tubulin acetylation inducer) is a highly potent and selective, reversible, cell-permeable inhibitor of HDAC6 ($IC_{50}=0.004\mu M$). IC_{50} 's for the other HDACs are 1000-fold higher, making tubacin both more selective and more potent than Tubastatin A, which also inhibits HDAC8. Concentration in cell culture experiments typically ranges from 2-50 μM .

Citations: 16

[View Online »](#)

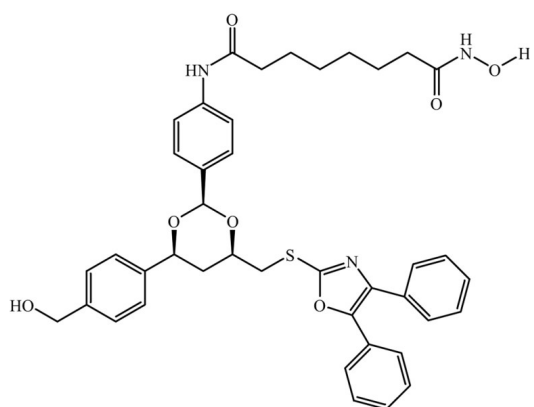
Ordering Information

[Order Online »](#)

BML-GR362-0500	500 μg
----------------	-------------

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 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	N-(4-((2R,4R,6S)-4-(((4,5-diphenyl-1,3-oxazol-2-yl)sulfanyl)methyl)-6-[4-(hydroxymethyl)phenyl]-1,3-dioxan-2-yl)phenyl)-N'-hydroxyoctanediamide
Appearance	White solid.
CAS	537049-40-4
Couple Target	HDAC, HDAC6
Couple Type	Inhibitor
Formula	C ₄₁ H ₄₃ N ₃ O ₇ S
MW	721.9
Purity	≥96% (ELSD)
Solubility	Soluble in DMSO (10 mg/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