

BML-281

HDAC6 inhibitor

Selective HDAC6 inhibitor ($IC_{50}=0.002nM$). IC_{50} values for other HDACs are 271, 252, 0.42, 6851 and 90.7 nM for HDACs 1,2,3,8 and 10 respectively. Potent inhibitor of the proliferation of a variety of pancreatic cancer cell lines (IC_{50} s in the 0.1-0.6 μ M range).

Citations: 2

[View Online »](#)

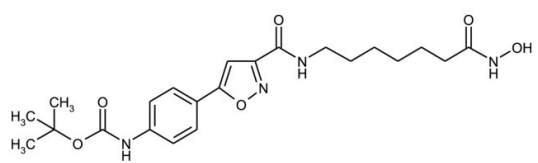
Ordering Information

[Order Online »](#)

BML-GR361-0001	1mg
BML-GR361-0005	5mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at -20°C for up to one year. Store solutions at -20°C for up to 3 months.
Short Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	N-Hydroxy-7-[5-(4-tertbutoxycarbonylaminophenyl)-3-isoxazolecarboxamido]heptanamide
Appearance	Off-white solid.
CAS	1045792-66-2
Couple Target	HDAC, HDAC6
Couple Type	Inhibitor
Formula	$C_{22}H_{30}N_4O_6$
MW	446.5
Purity	≥98% (TLC)
Solubility	Soluble in DMSO (15 mg/ml).

Last modified: May 29, 2024



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)