

MS-275

HDAC1 inhibitor

Preferentially inhibits HDAC1 (IC_{50} =300nM) over HDAC3 (IC_{50} =8 μ M). Has no inhibitory activity towards HDAC8 (IC_{50} >100 μ M).

Citations: 20

[View Online »](#)

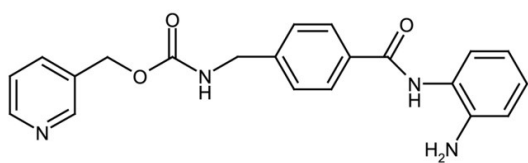
Ordering Information

[Order Online »](#)

ALX-270-378-M001	1mg
ALX-270-378-M025	25mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light. Packaged under inert gas.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	N-(2-Aminophenyl)-4-[N-(pyridine-3-ylmethoxy-carbonyl)aminomethyl]benzamide, Entinostat
Appearance	Off-white to tan solid.
CAS	209783-80-2
Couple Target	HDAC, HDAC1
Couple Type	Inhibitor
Formula	$C_{21}H_{20}N_4O_3$
MW	376.4
Purity	≥95% (1H-NMR)
Solubility	Soluble in DMSO 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