

Trichostatin C

HDAC inhibitor

Analog of trichostatin A (Prod. No. [BML-GR309](#)). Induces acetylation of histone H4. Shows antifungal activity.

Citations: 3

[View Online »](#)

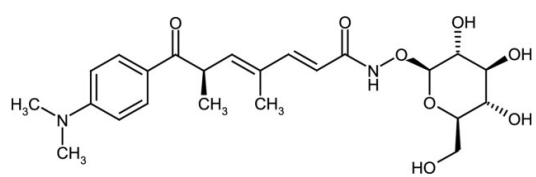
Ordering Information

[Order Online »](#)

ALX-380-239-MC05	0.5mg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Antibiotic 145-A

Appearance White to off-white powder.

CAS 68676-88-0

Couple Target HAT

Couple Type Activator

Formula $C_{23}H_{32}N_2O_8$

MI 14: 9649

MW 464.5

Purity ≥95% (HPLC)

Solubility Soluble in ethanol, methanol, DMF or DMSO. Limited water solubility.

Source Isolated from *Streptomyces* sp. MST-AS5376.

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-
be@enzolifesciences.com](mailto:info-be@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)