

Romidepsin

HDAC inhibitor

A potent and selective inhibitor of class I histone deacetylases (HDACs). Induces apoptosis in a variety of cell lines and displays antitumor activity in mouse models. Recently approved for clinical use in T-cell lymphoma.

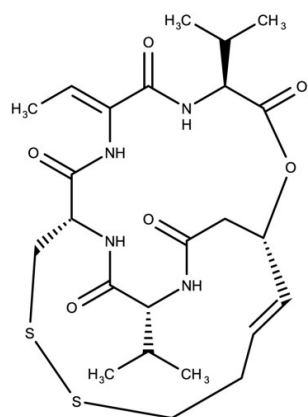
Ordering Information

[Order Online »](#)

ENZ-CHM468-0005	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 FK-228, FR901228, Depsipeptide

Appearance Colorless waxy solid – vials may appear empty to visual inspection.

CAS 128517-07-7

Couple Target HDAC

Couple Type Inhibitor

Formula $C_{24}H_{36}N_4O_6S_2$

Identity Determined by NMR.

MW 540.71

Purity ≥98% (HPLC)

Solubility Soluble in DMSO (10 mg/ml).

Last modified: February 27, 2025



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)