

Vorinostat

HDAC inhibitor, RNA splicing inhibitor.

Ordering Information

[Order Online »](#)

LKT-V5734-M100	100mg
LKT-V5734-M250	250mg
LKT-V5734-G001	1g

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Suberoylanilide hydroxamic acid, SAHA, Zolinza

CAS 149647-78-9

Couple Type Inhibitor

Formula $C_{14}H_{20}N_2O_3$

MW 264.32

Purity $\geq 98\%$



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