

Ebelactone A

Esterase inhibitor

Inhibitor of esterases, lipases and N-formylmethionine aminopeptidase.

Citations: 5

[View Online »](#)

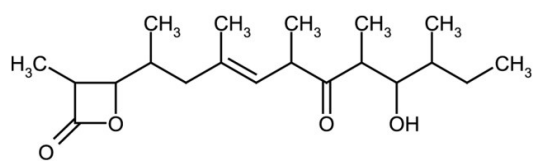
Ordering Information

[Order Online »](#)

ALX-270-013-M001	1mg
------------------	-----

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 3,11-Dihydroxy-2,4,6,8,10,12-hexamethyl-9-oxo-6-tetradecenoic 1,3-lactone

Appearance White crystalline powder.

CAS 76808-16-7

Couple Target Aminopeptidase, Esterase

Couple Type Inhibitor

Formula $C_{20}H_{34}O_4$

MW 338.5

Purity ≥99%

Solubility Soluble in methanol.

Source Microbial.

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