

Parthenolide

HDAC inhibitor

Sesquiterpene lactone. Widely used as an herbal remedy for arthritis and migraine. Has anti-inflammatory, antisecretory and spasmolytic activity. Specifically inhibits activation of NF- κ B by preventing the degradation of I κ B α and I κ B β . Inhibits activation of MAP kinase (MAPK/ERK) and generation of leukotriene B₄ and thromboxane B₂. Potent anticancer agent. Induces apoptosis in various cancer cell lines. Specifically inhibits histone deacetylase 1 (HDAC1) without affecting other class I/II HDACs. Inhibits tubulin carboxypeptidase (TCP) activity.

Citations: 22

[View Online »](#)

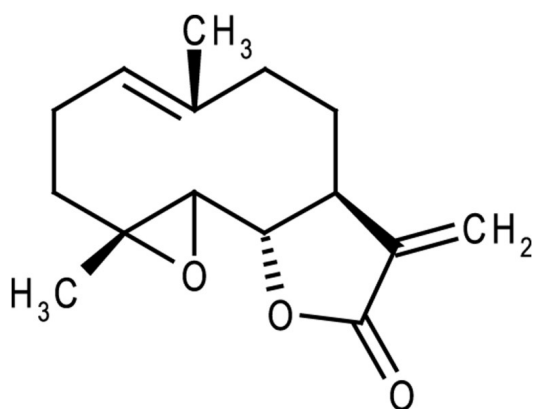
Ordering Information

[Order Online »](#)

BML-T113-0050	50mg
BML-T113-0250	250mg

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 (1aR-[1aR*,4E,7aS*,-10aS*,-10bR*])-2,3-6,7,7a,8,10a,10b-Octahydro-1a,5-dimethyl-8-methyleneoxireno[9,10]cyclo deca[1,2-b]furan-9(1aH)-one

Appearance White to off-white crystalline solid.

CAS 20554-84-1

Couple Target HDAC, HDAC1

Couple Type Inhibitor

Formula $C_{15}H_{20}O_3$

MI 14: 7048

MW 248.3

Purity ≥97% (HPLC, TLC)

RTECS LY4220000

Solubility Soluble in DMSO (100mg/ml), dichloromethane (50mg/ml) or 100% ethanol (20mg/ml).

Source Isolated from feverfew (*Chrysanthemum parthenium*) leaves.



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