

Embelin

XIAP inhibitor

Embelin is a novel cell-permeable inhibitor of XIAP (X-linked inhibitor of apoptosis) which was discovered by screening a library of natural products derived from traditional Chinese medicine. It binds to the XIAP BIR3 domain preventing its binding to and inhibition of caspase-9 and effector caspases. Embelin selectively inhibits cell growth in cancer cells expressing high levels of XIAP (IC₅₀ = 3-6 μ M) versus normal cells with low levels of XIAP.

Citations: 3

[View Online »](#)

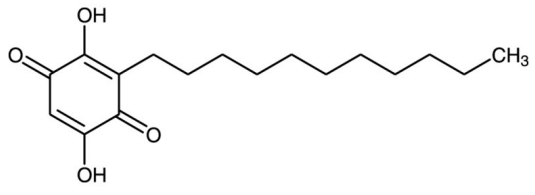
Ordering Information

[Order Online »](#)

BML-CM125-0010	10mg
----------------	------

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	Embelic acid
Appearance	Orange crystalline solid.
CAS	550-24-3
Couple Target	XIAP
Couple Type	Inhibitor
Formula	$C_{17}H_{26}O_4$
MW	294.4
Purity	≥98% (HPLC)
Solubility	Soluble in DMSO (10mg/ml)

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)