

Shikonin

Inhibits angiogenesis

Shikonin inhibits TNF- α - and B6 melanoma-induced angiogenesis in mice. It blocks expression of integrin $\alpha\beta 3$ and inhibits endothelial cell proliferation and migration *in vitro*.

Citations: 8

[View Online »](#)

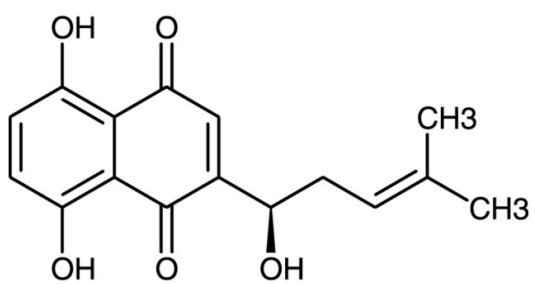
Ordering Information

[Order Online »](#)

BML-CT115-0010	10mg
BML-CT115-0050	50mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage +4°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance Red crystalline solid.

CAS 517-89-5

Couple Target CD51, CD61, Integrin

Couple Type Inhibitor

Formula $C_{16}H_{16}O_5$

Identity Determined by NMR and MS.

MW 288.3

Purity ≥98% (HPLC)

Solubility Soluble in DMSO.

Source Shikonin is a naphthoquinone isolated from the roots of *Lithospermum erythrorhizon*; a component of the traditional Chinese herbal remedy Shiunko.

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-
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)