

Wedelolactone

IKK inhibitor

Wedelolactone is a naturally occurring furanocoumarin isolated from *Eclipta prostrata*, an herbal medicine used to treat septic shock in traditional Chinese medicine. It inhibits the I κ B kinase (IKK) complex and suppresses LPS-induced caspase-11 expression in LPS-treated BALB 3T3 cells (IC₅₀ = 35 μ M)

Citations: 3

[View Online »](#)

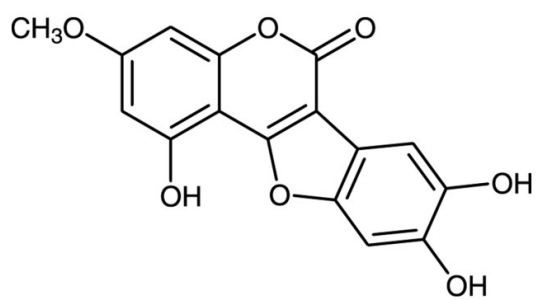
Ordering Information

[Order Online »](#)

BML-EI316-0001	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 7-Methoxy-5,11,12-trihydroxycoumestan

Appearance Yellow-green solid

CAS 524-12-9

Couple Target IKK, NF-kappaB

Couple Type Inhibitor

Formula $C_{16}H_{10}O_7$

MW 314.2

Purity ≥96% (HPLC)

Solubility Soluble in DMSO (25mg/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-
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)