

Tanshinone IIA

AP-1 inhibitor

Tanshinone IIA is a diterpenoid naphthoquinone found in the traditional Chinese medicine Tanshen (*Salvia sp*). It inhibits AP-1 activity by suppressing jun-fos-DNA complex formation ($IC_{50}=0.22\text{ }\mu\text{M}$) and induces apoptosis in a variety of cell lines

Citations: 4

[View Online »](#)

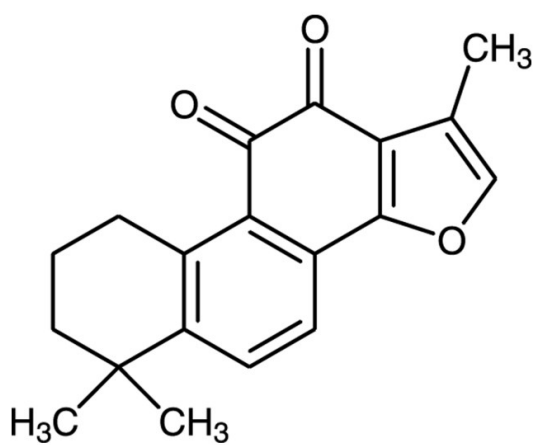
Ordering Information

[Order Online »](#)

BML-GR336-0025	25mg
----------------	------

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 Orange solid.

CAS 568-72-9

Couple Target Jun

Couple Type Inhibitor

Formula $C_{19}H_{18}O_3$

MW 294.3

Purity ≥93% (HPLC)

Solubility Soluble in DMSO (25mg/ml) or 100% ethanol (5mg/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)