

Hinokitiol

Chelator that can induce differentiation in tumor cells

Hinokitiol (β -thujaplicin) is a potent metal chelator that induces differentiation and apoptosis in teratocarcinoma F9 cells and tyrosinase, most likely *via* metal chelation.

Citations: 6

[View Online »](#)

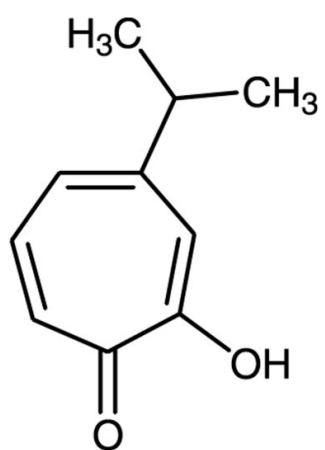
Ordering Information

[Order Online »](#)

BML-FR110-0250	250mg
----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

CAS	499-44-5
MW	164.2
Purity	≥98% (GC)

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)