

Ellipticine

Topoisomerase inhibitor

A DNA-intercalating antitumor agent. Inhibits DNA topoisomerase II. Inhibits the phosphorylation of p53 in Lewis lung carcinoma cells and induces apoptosis. Inhibits cytochrome P-450-1A1.

Citations: 4

[View Online »](#)

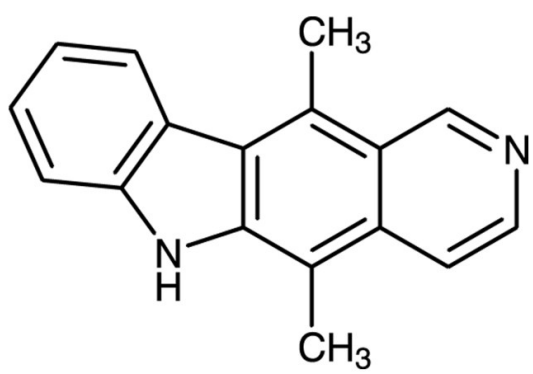
Ordering Information

[Order Online »](#)

BML-GR315-0010	10mg
BML-GR315-0050	50mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 5,11-Dimethyl-6H-pyrido[4,3-b]carbazole

Appearance Bright yellow solid.

CAS 519-23-3

Couple Target Topoisomerase

Couple Type Inhibitor

Formula $C_{17}H_{14}N_2$

Identity Determined by NMR.

MW 246.3

Purity ≥98% (TLC)

Solubility Soluble in DMSO, DMF or ethanol.

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)