

3,3'-Diindolylmethane

Cytochrome P450 activator

Dimer of indole-3-carbinol (Prod. No. ALX-350-347). Anticancer and antineoplastic agent. Induces apoptosis in human cancer cells. Activator of cytochrome P450. Partial agonist of aryl hydrocarbon receptor (AhR).

Citations: 10

[View Online »](#)

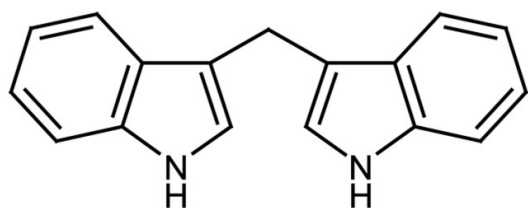
Ordering Information

[Order Online »](#)

ALX-350-370-M500	500mg
------------------	-------

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 DIM

Appearance White to light pink powder.

CAS 1968-05-4

Couple Target Cytochrome P450

Couple Type Activator

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

MW 246.3

Purity ≥98% (HPLC)

Solubility Soluble in DMSO (200mg/ml); insoluble in water.

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)