

2-Methoxyestradiol

Estrogen receptor inhibitor

An endogenous estrogen metabolite with low affinity for the estrogen receptor but which disrupts microtubule function. Potent inhibitor of endothelial cell proliferation and migration. Inhibits bFGF and VEGF-induced corneal neovascularization in mice. Posttranscriptionally downregulates HIF-1 α expression.

Citations: 3

[View Online »](#)

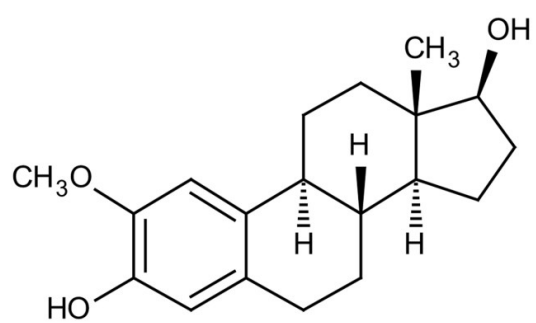
Ordering Information

[Order Online »](#)

BML-S540-0010	10mg
BML-S540-0050	50mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 2-ME

Appearance White solid.

CAS 362-07-2

Couple Target Estrogen receptor

Couple Type Inhibitor

Formula $C_{19}H_{26}O_3$

Identity Determined by ¹H-NMR and MS.

MW 302.4

Purity ≥95% (TLC)

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)