

Gossypol

Male antifertility agent

Mixture of ~65% (+)- and ~35% (-)-enantiomers. Male antifertility agent with antispermatogenic activity. Shows antitumor, antiviral and antioxidant actions. Small molecule inhibitor of Bcl-2/Bcl-X_L. Induces apoptosis. Synergizes with radiation and chemotherapy *in vitro*. Reversible inhibitor of protein phosphatase 2B (PP2B; calcineurin). Protein kinase C (PKC) inhibitor. Inhibits activities of the mitotic kinesin Eg5 (IC₅₀=10.8μM/L). Inhibits RBL-1, 5-lipoxygenase (IC₅₀=0.3μM) and 12-lipoxygenase (IC₅₀=0.7μM). Inhibits PAF and leukotriene-induced guinea pig parenchyma contractions. Novel direct-acting inhibitor of calcineurin.

Citations: 21

[View Online »](#)

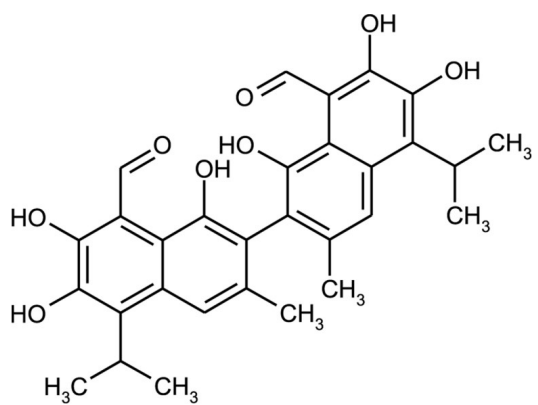
Ordering Information

[Order Online »](#)

BML-EI130-0100	100mg
----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored at -20°C. Stock solutions are stable for up to 1 month at -20°C. Avoid alkaline solutions.
Handling	Protect from light and oxygen. Keep under inert gas.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	2,2'-bis[8-Formyl-1,6,7-trihydroxy-5-isopropyl-3-methylnaphthalene]
Appearance	Off-white to yellow solid.
CAS	303-45-7
Couple Target	Kinesin, Lipoygenase, PKC, Serine/threonine-protein phosphatase
Couple Type	Inhibitor
Formula	$C_{30}H_{30}O_8$
MI	14: 4528
MW	518.6
Purity	≥90% (HPLC)
RTECS	DU3100000
Solubility	Soluble in 100% ethanol (25mg/ml), dimethylformamide (25mg/ml), acetone or DMSO; sparingly soluble in methanol (2mg/ml); insoluble in water.

Source Isolated from *Gossypium* genus, Malvaceae.



Technical Info / Product

ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +49 7621 5500 526
info@enzolifesciences.com
eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
fr@enzolifesciences.com
be@enzolifesciences.com

France
Phone: +33 472 440 655
info@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info@enzolifesciences.com
de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info@enzolifesciences.com
uk@enzolifesciences.com

Notes

Note: Product is not sterile.