

# 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<sub>L</sub>. 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 (IC<sub>50</sub>=10.8μM/L). Inhibits RBL-1, 5-lipoxygenase (IC<sub>50</sub>=0.3μM) and 12-lipoxygenase (IC<sub>50</sub>=0.7μM). Inhibits PAF and leukotriene-induced guinea pig parenchyma contractions. Novel direct-acting inhibitor of calcineurin.

Citations: 21

[View Online »](#)

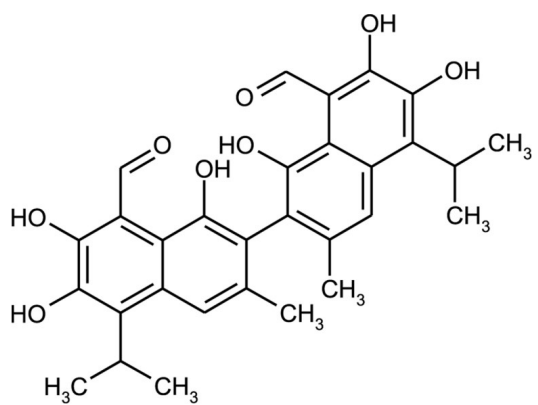
## Ordering Information

[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, Lipoxygenase, PKC, Serine/threonine-protein phosphatase
Couple Type	Inhibitor
Formula	C <sub>30</sub> H <sub>30</sub> O <sub>8</sub>
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 <i>Gossypium</i> genus, Malvaceae.

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)