

# Gambogic acid

Bcl-2 inhibitor, Bax activator

Gambogic Acid is a natural product isolated from *Garcinia Hurburyi*. It induces apoptosis in several tumor cell lines including T47D breast cancer cells and MGC-803 human gastric carcinoma cells but not in normal cells. Apoptosis induction proceeds *via* upregulation of bax and inhibition of bcl-2 expression. The target of gambogic acid has been identified as the transferrin receptor.

Citations: 3

[View Online »](#)

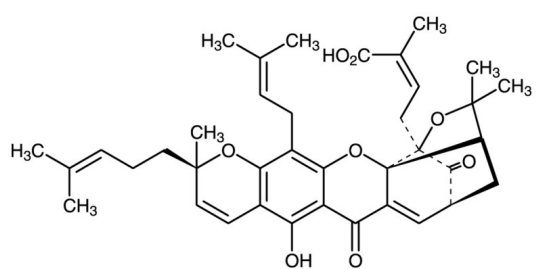
## Ordering Information

[Order Online »](#)

BML-AP305-0005	5mg
BML-AP305-0025	25mg

## 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

**Appearance** Yellow-orange solid

**CAS** 2752-65-0

**Couple Target** Bax, Bcl-2

**Couple Type** Activator, Inhibitor

**Formula**  $C_{38}H_{44}O_8$

**MW** 628.8

**Purity** ≥95% (TLC)

**Solubility** Soluble in DMSO (25 mg/ml) or ethanol (25 mg/ml)

**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)