

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 »](#)

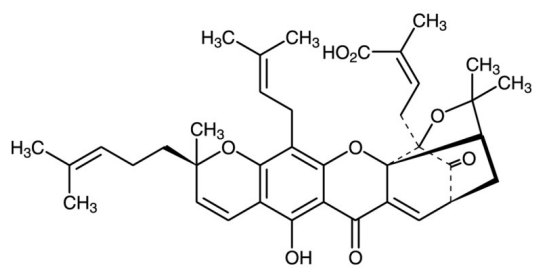
Ordering Information

[Order Online »](#)

BML-AP305-0005	5mg
----------------	-----

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

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

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

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

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

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