

Garcinol

HAT inhibitor

Garcinol is a polyisoprenylated benzophenone derivative isolated from *Garcinia indica*. It is a potent inhibitor of histone acetyltransferases (HATs) p300 ($IC_{50}=7\mu M$) and PCAF ($IC_{50}=5\mu M$) both *in vitro* and *in vivo*. HAT activity-dependent chromatin transcription was inhibited by garcinol, but it had no effect on naked DNA transcription. It induces apoptosis in HeLa cells ($100\mu M$) where it down regulates global gene expression. Acetylation of histone H4 by p300 was more sensitive to inhibition by garcinol compared with that of H3. Histone H4 acetylation was completely inhibited at $1\mu M$ while acetylation of H3 was not abolished at $20\mu M$.

Citations: 18

[View Online »](#)

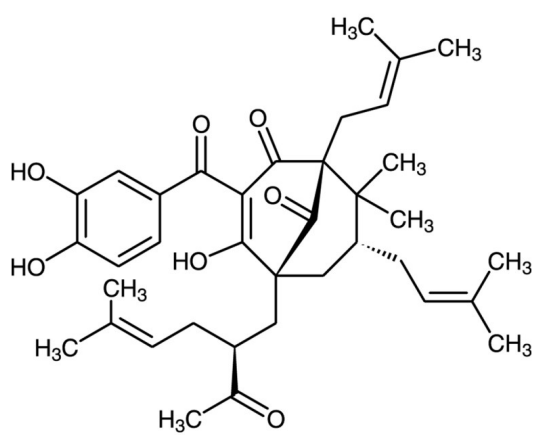
Ordering Information

[Order Online »](#)

BML-GR343-0010	10mg
BML-GR343-0050	50mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at -20°C for up to 1 year. Store solutions at -20°C for up to 3 months.
Handling	Protect from exposure to air
Short Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	Yellow solid.
CAS	78824-30-3
Couple Target	HAT
Couple Type	Inhibitor
Formula	$C_{38}H_{50}O_6$
Identity	Determined by NMR.
MW	602.8
Purity	≥95% (TLC)
Solubility	Soluble in DMSO (25mg/ml) or 100% ethanol (25mg/ml) or dimethylformamide (25mg/ml).
Source	Synthetic.

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-
be@enzolifesciences.com](mailto:info-be@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)