

Geranylgeraniol

Geranylgeranylpyrophosphate precursor;
apoptosis inducer

Precursor to geranylgeranylpyrophosphate. Induces apoptosis in HL-60 cells.

Citations: 1

[View Online »](#)

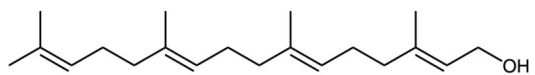
Ordering Information

[Order Online »](#)

BML-G243-0020	20mg
---------------	------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Keep under inert gas. Light and Temperature sensitive.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Appearance	Colorless Oil.
CAS	24034-73-9
Formula	$C_{20}H_{34}O$
Identity	Determined by 1H-NMR.
MW	290.5
Purity	≥95% (TLC)
Solubility	Soluble in chloroform or ethyl acetate.

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)