

Gingerol

ATPase activator

Gingerol ([6]-gingerol), a pungent vanilloid isolated from ginger rhizomes (*Zingiber officinale*), is a novel and selective activator of skeletal and cardiac SR Ca²⁺-ATPase (EC₅₀=4μM). It is a novel cardiogenic agent with antiplatelet activity. It induces apoptosis in HL-60 cells. Acts as an agonist at vanilloid receptors.

Citations: 6

[View Online »](#)

Ordering Information

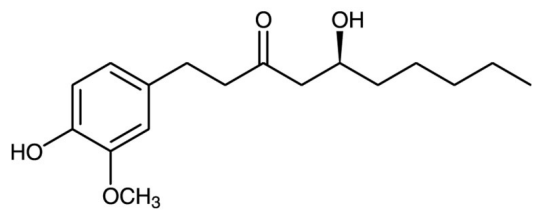
[Order Online »](#)

BML-CA422-0005

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	White to pale yellow oily or waxy solid.
CAS	23513-14-6
Couple Target	ATPase
Couple Type	Activator
Formula	$C_{17}H_{26}O_4$
MW	294.4
MeltingPoint	31°C
Purity	≥97% (HPLC)
Solubility	Soluble in 100% ethanol.

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)