

Ilimaquinone

ADP-ribosylation factor inhibitor

Displays antimicrobial, anti-HIV and anti-inflammatory properties. Induces complete and reversible breakdown of Golgi membranes. Inhibits the association of β -COP and ADP-ribosylation factor to Golgi membranes, but unlike BFA, does not induce retrograde transport of Golgi enzymes into ER or cause random fusion of successive Golgi cisternae. Interacts with methylation enzymes. Shows anticancer activity.

Citations: 8

[View Online »](#)

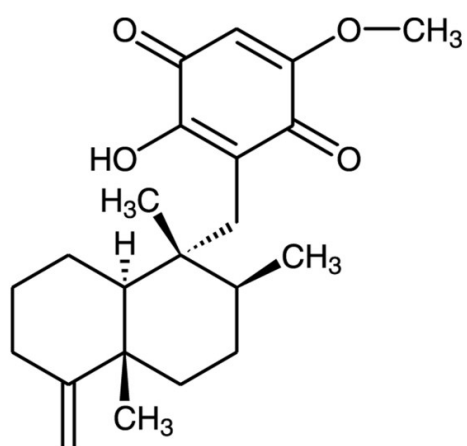
Ordering Information

[Order Online »](#)

BML-G435-0100	100 μ g
---------------	-------------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Solutions can be stored at -20°C for up to 3 months.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	3-[(Decahydro-1β,2β,4αβ-trimethyl-5-methylene-1-naphthyl)methyl]-2-hydroxy-5-methoxybenzoquinone
Appearance	Yellow oil.
CAS	71678-03-0
Couple Target	ADP-ribosylation factor
Couple Type	Inhibitor
Formula	$C_{22}H_{30}O_4$
MW	358.5
Purity	≥98% (TLC)
Solubility	25mg/ml soluble in DMSO or 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)