

Ginkgolide B

PAF receptor ligand

Most potent member of the ginkgolide family of PAF antagonists. Ginkgolide B accelerates PAF hydrolysis by promoting PAF-AH I α_2 homodimer activity. Inhibits PAF-induced PMNL chemotaxis (IC_{50} =490 nM). Orally effective against PAF-induced bronchospasm.

Citations: 6

[View Online »](#)

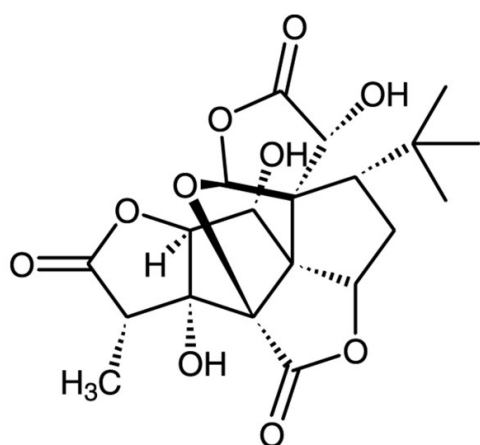
Ordering Information

[Order Online »](#)

BML-L135-0010	10mg
---------------	------

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

Alternative Name BN 52021

Appearance White crystalline powder.

CAS 15291-77-7

Couple Target PAF receptor

Couple Type Ligand

Formula $C_{20}H_{24}O_{10}$

MW 424.4

Purity ≥90% (HPLC)

Solubility Soluble in DMSO (25mg/ml) or 100% ethanol (2mg/ml).

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-
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)