

Brefeldin A

GEF inhibitor

Inhibitor of intracellular protein transport and protein secretion, interfering with trafficking in the *trans*-Golgi network, leading to the accumulation of cycling proteins in ERGIC clusters. Blocks ADP-ribosylation factor (Arf) in an inactive GDP-bound conformation and thereby prevents binding of COPI coats to ERGIC and Golgi membranes. Upon brefeldin A treatment the Golgi rapidly tubulates and fuses with the ER by an energy-, temperature-, and microtubule-dependent process. In contrast, the drug has little effect on the ERGIC, which keeps its identity, although the ERGIC clusters become larger and more uniformly distributed in the cytoplasm of the cells. Inhibits intracellular collagen degradation. Also inhibits apical Na⁺ channels in epithelia. Induces apoptosis.

Citations: 39

[View Online »](#)

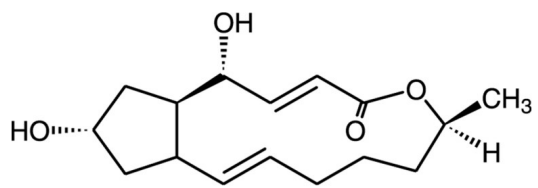
Ordering Information

[Order Online »](#)

BML-G405-0005	5mg
BML-G405-0025	25mg

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 BFA, Ascotoxin, Decumbin

Appearance White to off-white solid.

CAS 20350-15-6

Couple Target GEF

Couple Type Inhibitor

Formula $C_{16}H_{24}O_4$

MI **14:** 1369

MW 280.4

Purity ≥98% (HPLC)

Solubility Soluble in DMSO (20mg/ml), 100% ethanol (5mg/ml), methanol, acetone or ethyl acetate (1mg/ml).

Source Isolated from *Eupenicillium brefeldianum*.

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)