

Fusicoccin

Plant H⁺-ATPase activator

Fusicoccin is a diterpene glucoside with numerous effects. Fusicoccin induces H⁺ extrusion from plant cells by 14-3-3-dependent activation of the plasma membrane H⁺-ATPase. Fusicoccin induces translocation of 14-3-3 proteins from the cytosol to the membrane and causes a dramatic increase in cytosolic 14-3-3. Fusicoccin stimulates release of cytochrome c from mitochondria, cell enlargement, ion uptake, seed germination, opening the stomata, etc. Fusicoccin markedly reduces the ability of the cells to scavenge H₂O₂. Fusicoccin has been demonstrated to cause randomization of asymmetry of the left-right axis (heterotaxia) in *Xenopus* embryos.

Citations: 16

[View Online »](#)

Ordering Information

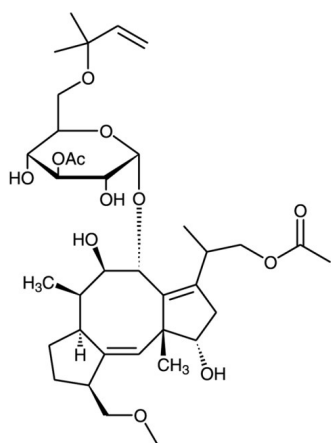
[Order Online »](#)

| | |
|----------------|-----|
| BML-EI334-0001 | 1mg |
|----------------|-----|

Manuals, SDS & CofA

[View Online »](#)

- Phytotoxin used in plant research



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 NSC 113500

Appearance Solid.

CAS 20108-30-9

Couple Target 14-3-3, ATPase

Couple Type Activator, Inhibitor

Formula $C_{36}H_{56}O_{12}$

MW 680.8

Purity ≥90% (HPLC)

Solubility Soluble in DMSO.



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)