

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: 17

[View Online »](#)

Ordering Information

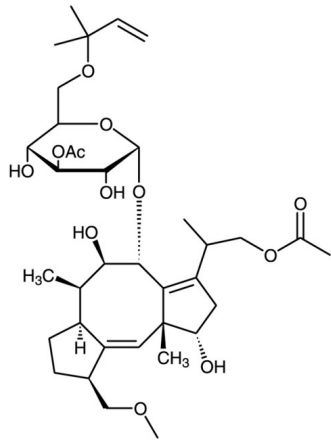
[Order Online »](#)

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	Film.
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.

Last modified: August 15, 2025



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com