

# Fusicoccin

## Plant H<sup>+</sup>-ATPase activator

Fusicoccin is a diterpene glucoside with numerous effects. Fusicoccin induces H<sup>+</sup> extrusion from plant cells by 14-3-3-dependent activation of the plasma membrane H<sup>+</sup>-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 H<sub>2</sub>O<sub>2</sub>. Fusicoccin has been demonstrated to cause randomization of asymmetry of the left-right axis (heterotaxia) in *Xenopus* embryos.

Citations: 17

[View Online »](#)

## Ordering Information

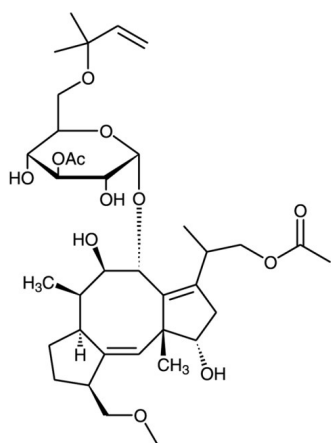
[Order Online »](#)

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](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)