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 »

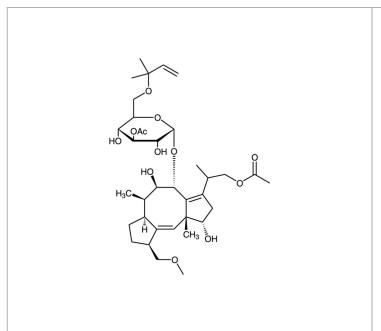
BML-EI334-0001

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.

Last modified: August 15, 2025

