

# Anisomycin

Activator of MAPK, Jnk and p38

Antibiotic. Activates JNK/SAPKs, p38 and MAP kinases but not ERKs leading to the phosphorylation of S6, histone H3 and HMG-14 and the induction of the immediate-early genes c-fos and c-jun. Induces apoptosis in the human monoblastoid cell line. Used at 300 ng/ml. Inhibits peptidyl transferase at higher concentrations. Used in the eradication of bean mildew. Inhibits other pathogenic fungi in plants.

Citations: 18

[View Online »](#)

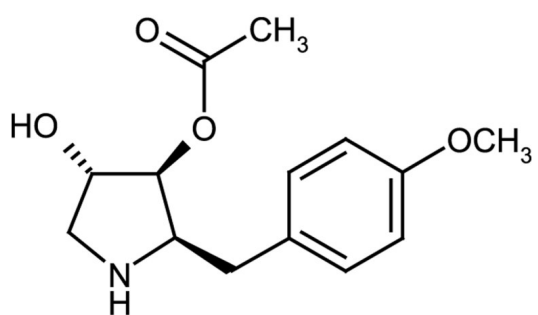
## Ordering Information

[Order Online »](#)

BML-ST102-0050	50mg
----------------	------

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored, as supplied, at -20°C. Stock solutions are stable for up to 3 months at -20°C.
Long Term Storage	-20°C
Shipping	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	2-(p-Methoxybenzyl)-3,4-pyrrolidinediol-3-acetate
Appearance	White solid.
CAS	22862-76-6
Couple Target	Jnk, MAPK, p38
Couple Type	Activator
Formula	$C_{14}H_{19}NO_4$
MI	14: 670
MW	265.3
Purity	≥97% (TLC)
RTECS	BZ9800000
Solubility	Soluble in DMSO (25mg/ml).
Source	Isolated from <i>Streptomyces griseolus</i> .



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)