

# Citrinin

## MPTP activator

Antibiotic. Induces mitochondrial permeability pore opening and inhibits respiration by interfering with complex I of the respiratory chain. Mycotoxin. Acts as nephrotoxin in all species in which it has been tested. Has been implicated as a cause of Balkan nephropathy and yellow rice fever in humans. Induces apoptosis.

Citations: 5

[View Online »](#)

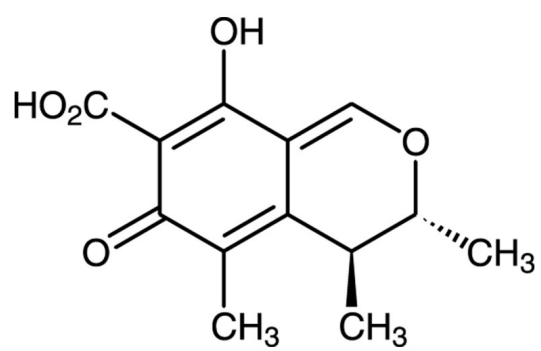
## Ordering Information

[Order Online »](#)

BML-CM116-0005	5mg
BML-CM116-0025	25mg

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Solutions change color with changes in pH from yellow (pH 4.6) to red (pH 9.9).
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

Appearance	Yellow crystalline solid.
CAS	518-75-2
Couple Target	MPTP
Couple Type	Activator
Formula	$C_{13}H_{14}O_5$
MW	250.2
Purity	≥98% (HPLC, TLC)
Solubility	Soluble in 100% ethanol, methylene chloride (10mg/ml), methanol, dioxane, pyridine or dichloromethane. Insoluble in water.
Source	Isolated from <i>Penicillium citrinum</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-  
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)