

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 »](#)

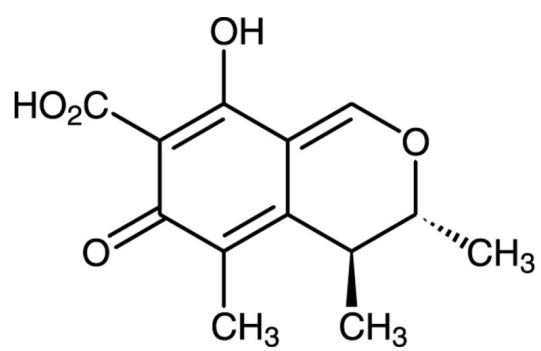
Ordering Information

[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 dichlormethane. Insoluble in water.
Source	Isolated from <i>Penicillium citrinum</i>

Last modified: May 29, 2024



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)