

Ochratoxin A

Mycotoxin. Nephrotoxic.

Mycotoxin. Natural contaminant of mouldy food and feed. It has a number of toxic effects, the most prominent being nephrotoxicity. Immunosuppressive, genotoxic, teratogenic and carcinogenic. Stimulates lipid peroxidation. Induces oxidative DNA lesions coupled with direct DNA adducts via quinone formation.

Citations: 13

[View Online »](#)

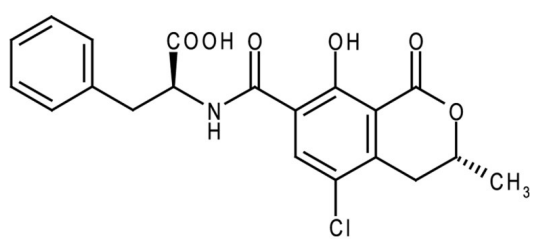
Ordering Information

[Order Online »](#)

ALX-630-089-M001	1mg
ALX-630-089-M005	5mg
ALX-630-089-M025	25mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Long Term Storage +4°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance White to off-white crystalline powder.

CAS 303-47-9

Formula $C_{20}H_{18}ClNO_6$

MI **14:** 6740

MW 403.8

Purity ≥98% (HPLC, TLC)

RTECS AY4375000

Solubility Soluble in methanol (10mg/ml).

Source Isolated from *Aspergillus ochraceus*.

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)