

Aflatoxin G2

Hepatotoxic and hepatocarcinogenic
mycotoxin

Potent hepatotoxic and hepatocarcinogenic mycotoxin.

Not for sale in Turkey and South Africa.

May require a license for import, please [contact us](#) for more information.

Citations: 3

[View Online »](#)

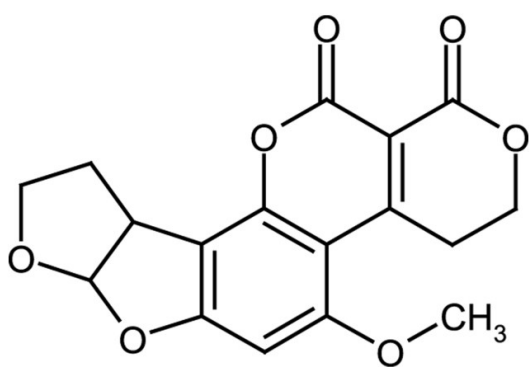
Ordering Information

[Order Online »](#)

ALX-630-106-M001	1mg
ALX-630-106-M005	5mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Handling Protect from light.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance White to yellow powder. Blue-green fluorescence.

CAS 7241-98-7

Formula $C_{17}H_{14}O_7$

MI **14:** 181

MW 330.3

Purity ≥98% (TLC)

RTECS LV1700000

Solubility Soluble in methylene chloride (5mg/ml), DMSO or methanol.

Source Isolated from *Aspergillus flavus*.

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-
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)