

Aflatoxin M₂

Hepatotoxic and hepatocarcinogenic mycotoxin

Potent hepatotoxic and hepatocarcinogenic mycotoxin. Metabolite of aflatoxin B₂. DNA-damaging agent.

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

Citations: 4

[View Online »](#)

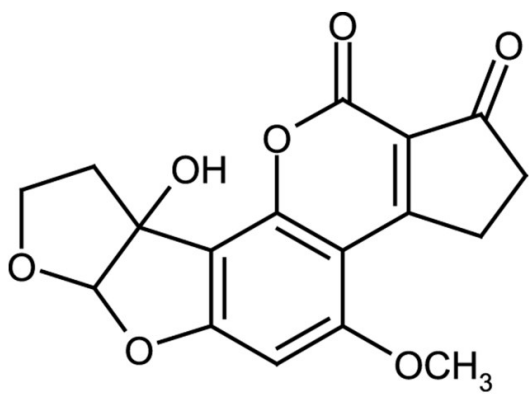
Ordering Information

[Order Online »](#)

ALX-630-114-MC01	0.1mg
------------------	-------

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

Alternative Name	4-Hydroxyafatoxin B2
Appearance	Off-white powder.
CAS	6885-57-0
Formula	$C_{17}H_{14}O_7$
MI	14: 182
MW	330.3
Purity	≥98% (HPLC, TLC)
RTECS	GY1720000
Solubility	Soluble in DMSO or methanol; insoluble in water.
Source	Isolated from <i>Aspergillus flavus</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)