

Aflatoxin B1

Mycotoxin

Naturally occurring mycotoxin produced by many species of *Aspergillus*. Metabolized by the liver to the reactive intermediate aflatoxin M₁. Exposure may produce acute necrosis, cirrhosis and carcinoma of the liver. Induces DNA damage.

Not for sale in Turkey and South Africa.

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

Citations: 10

[View Online »](#)

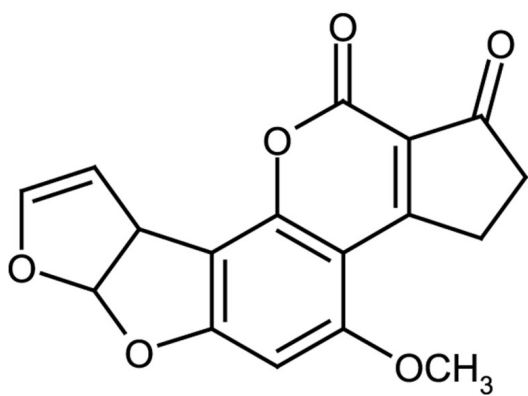
Ordering Information

[Order Online »](#)

ALX-630-093-M001	1mg
ALX-630-093-M005	5mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. May be stored in solution for at least three to four weeks at -20°C.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	White to yellow powder.
CAS	1162-65-8
Formula	$C_{17}H_{12}O_6$
MI	14: 180
MW	312.3
Purity	≥98% (HPLC, TLC)
RTECS	GY1925000
Solubility	Soluble in DMSO (20mg/ml) or methylene chloride (10mg/ml).
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)