

Aflatoxin M₁

Hepatotoxic and hepatocarcinogenic mycotoxin

Potent hepatotoxic and hepatocarcinogenic mycotoxin. Hydroxylated metabolite of aflatoxin B₁. Maximum absorbance (λ_{max}) at 228, 264 and 360nm. *Not for sale in Turkey and South Africa. May require a license for import, please [contact us](#) for more information.*

Citations: 4

[View Online »](#)

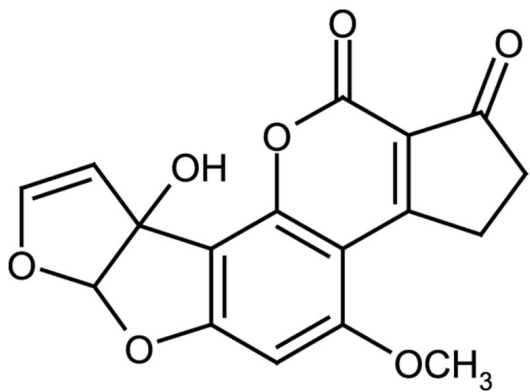
Ordering Information

[Order Online »](#)

ALX-630-095-MC01	0.1mg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance White to yellow powder.

CAS 6795-23-9

Formula $C_{17}H_{12}O_7$

MI **14:** 182

MW 328.3

Purity ≥95% (HPLC, TLC)

RTECS GY1880000

Solubility Soluble in methylene chloride (1mg/ml).

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)