

# Aflatoxin M1

Hepatotoxic and hepatocarcinogenic mycotoxin

Potent hepatotoxic and hepatocarcinogenic mycotoxin. Hydroxylated metabolite of aflatoxin B<sub>1</sub>. Maximum absorbance ( $\lambda_{\text{max}}$ ) 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 »](#)

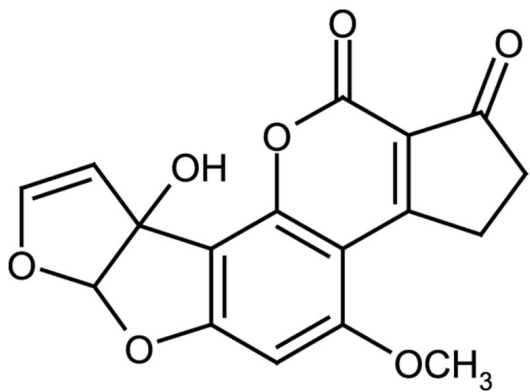
Ordering Information

[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	+4°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	<b>14:</b> 182
MW	328.3
Purity	≥95% (HPLC, TLC)
RTECS	GY1880000
Solubility	Soluble in methylene chloride (1mg/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-  
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)