

Sterigmatocystin

Mycotoxin

Mycotoxin produced by strains of the common molds. Inhibitor of DNA synthesis. Induces sister chromatid exchanges in bone marrow cells of mice. Causes necrosis.

Citations: 7

[View Online »](#)

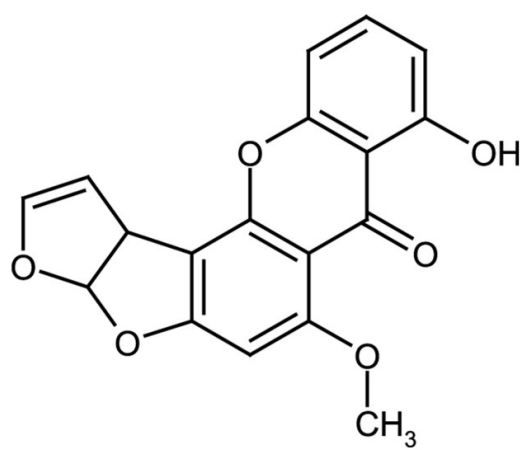
Ordering Information

[Order Online »](#)

ALX-630-116-M001	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	3a,12c-Dihydro-8-hydroxy-6-methoxy-7H-furo[3',2':4,5]furo[2,3-c]xanthen-7-one
Appearance	Pale yellow solid.
CAS	10048-13-2
Formula	$C_{18}H_{12}O_6$
MW	324.3
Purity	≥98% (TLC, HPLC)
RTECS	LV1750000
Solubility	Soluble in DMSO (10mg/ml), chloroform (71mg/ml), pyridine (18mg/ml) or acetone (5mg/ml).
Source	Isolated from <i>Aspergillus versicolor</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)