

Methoxsalen

Cytochrome P450 inhibitor

A potent suicide inhibitor of cytochrome P450. Antipsoriatic activity following ultraviolet light exposure.

Citations: 3

[View Online »](#)

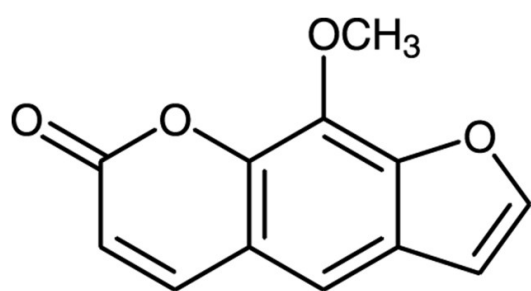
Ordering Information

[Order Online »](#)

BML-EI132-0001	1g
----------------	----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Xanthotoxin

Appearance White to yellow powder.

CAS 298-81-7

Couple Target Cytochrome P450

Couple Type Inhibitor

Formula $C_{12}H_8O_4$

MW 216.2

Purity $\geq 98\%$ (GC)

Solubility Soluble in acetone or hot ethanol (slight).

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)