

Triacsin C

Acyl-CoA synthetase inhibitor

Analog of polyunsaturated fatty acid. Potent inhibitor of long-chain fatty acyl CoA synthetase. Selectively inhibits arachidonoyl-CoA synthetase in intact cells and the nonspecific acyl-CoA synthetase in cell sonicates $IC_{50}=3.6-8.7\mu M$. Inhibits neutrophil functions. Blocks β -cell apoptosis induced by fatty acids (lipoapoptosis) and synthesis of triglycerides, diglycerides and cholesterol esters. Potent vasodilator.

Citations: 22

[View Online »](#)

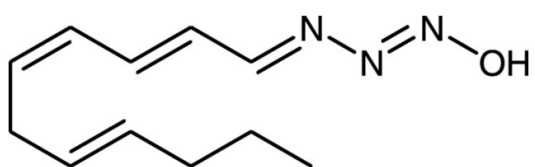
Ordering Information

[Order Online »](#)

BML-EI218-1000	1mg
BML-EI218-0100	100 μ g

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for up to 1 year after receipt when stored at -20°C.
Handling	Protect from light. Keep under inert gas.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	2,4,7-Undecatrienal nitrosohydrazone, 2E,4E,7E-Undecatriene-1-triazene, WS1228A
Appearance	Liquid. In DMSO.
CAS	76896-80-5
Couple Target	Acyl-CoA synthetase
Couple Type	Inhibitor
Formula	$C_{11}H_{17}N_3O$
MW	207.3
Purity	≥90% (TLC)
Solubility	Soluble in DMSO (25mg/ml) or methanol (5mg/ml).
Technical Info / Product Notes	Note: Product is not sterile.

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)