

Cerulenin

Fatty acid synthase inhibitor

Antibiotic. Fatty acid synthase (FAS) inhibitor reported to bind in equimolar ratio to β -keto-acyl-ACP synthase (KAS; $IC_{50}=1.5\mu M$), thus inhibiting protein acylation at concentrations of 45-134 μM . Produces metabolic effects similar to the effects of leptin, but through mechanisms that are independent of, or down-stream from, both leptin and melanocortin receptors. Apoptosis inducer. Inhibitor of bacterial fatty acid synthesis (inhibits FabH, FabB and FabF condensation enzymes).

Citations: 11

[View Online »](#)

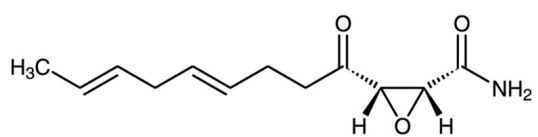
Ordering Information

[Order Online »](#)

BML-G237-0005	5mg
BML-G237-0025	25mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored, as supplied, at -20°C. Stock solutions are stable for up to 3 months at -20°C.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	2,3-Epoxy-4-oxo-7,10-dodecadienamide
Appearance	White to off-white solid.
CAS	17397-89-6
Couple Target	Fatty acid synthase
Couple Type	Inhibitor
Formula	$C_{12}H_{17}NO_3$
MI	14: 2004
MW	223.3
Purity	≥98% (TLC)
RTECS	JR1670000
Solubility	Soluble in DCM (5mg/ml).



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)