

Melittin (natural)

PLA₂ activator

Causes disruption of normal cellular activity and cell lysis. Binds calmodulin in a calcium-dependent manner. Activates phospholipase A₂ (PLA₂) and inhibits protein kinase C (PKC) by binding to the catalytic domain in a Mg-ATP sensitive manner. Antirheumatic.

Citations: 10

[View Online »](#)

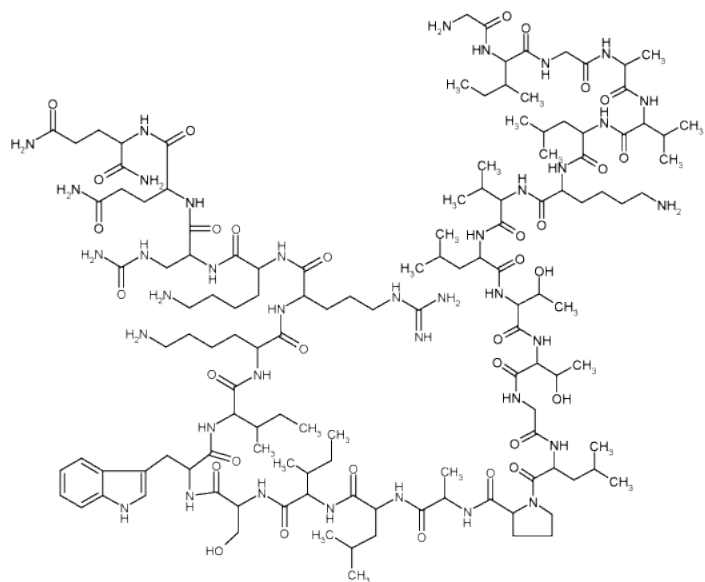
Ordering Information

[Order Online »](#)

ALX-162-006-M005	5mg
------------------	-----

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Very hygroscopic. Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Appearance	White to faint tan powder.
CAS	20449-79-0
Formula	$C_{131}H_{229}N_{39}O_{31}$
MI	14: 5824
MW	2846.5
Purity	≥85% (HPLC)
Purity Detail	Phospholipase A2 impurity:
RTECS	OS3960000
Sequence	H-Gly-Ile-Gly-Ala-Val-Leu-Lys-Val-Leu-Thr-Thr-Gly-Leu-Pro-Ala-Leu-Ile-Ser-Trp-Ile-Lys-Arg-Lys-Arg-Gln-Gln-NH ₂
Solubility	Soluble in PBS (5mg/ml) or water (5mg/ml).
Source	Isolated from bee venom.

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-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@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)