

Melittin (synthetic)

PLA₂ activator

PLA₂ stimulatory bee venom peptide. Sequence is similar to Phospholipase A₂ Activating Peptide (Prod. No. ALX-165-007). Activates PLA₂ (5-fold at 10µg/ml) in BC3H1 cells. For additional technical information see melittin (natural) (Prod. No. ALX-162-006).

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-162-007-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

Appearance	White to off-white powder.
CAS	20449-79-0
Formula	$C_{131}H_{229}N_{39}O_{31}$
MI	14: 5824
MW	2846.5
Peptide Content	70-90%
Purity	≥97% (HPLC)
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 water.
Source	Synthetic.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com