

Gramicidin A

Ionophore antibiotic

Naturally occurring ion channel forming pentadecapeptide. Causes K^+/H^+ -exchange in mitochondria in a non-voltage dependent manner.

Citations: 6

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-350-233-M001	1mg
ALX-350-233-M005	5mg

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

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

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance White to off-white powder.

CAS 11029-61-1

Formula $C_{99}H_{140}N_{20}O_{17}$

MI **14:** 4532

MW 1882.3

Purity ≥90% (HPLC)

Purity Detail Contaminants: ≤5% Gramicidin C (HPLC).

RTECS MD8230000

Sequence HCO-Val-Gly-Ala-D-Leu-Ala-D-Val-Val-D-Val-Trp-D-Leu-Trp-D-Leu-Trp-D-Leu-Trp-NHCH₂CH₂OH

Solubility Soluble in DMSO, lower alcohols or acetic acid.

Source Isolated from *Bacillus brevis*.

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)