

Cyclo [Arg-Ala-Asp-D-Phe-Val]

Negative control for Cyclo [Arg-Gly-Asp-D-Phe-Val]

Negative control for integrin $\alpha\text{v}\beta 3$ antagonist Cyclo [Arg-Gly-Asp-D-Phe-Val].

Citations: 1

[View Online »](#)

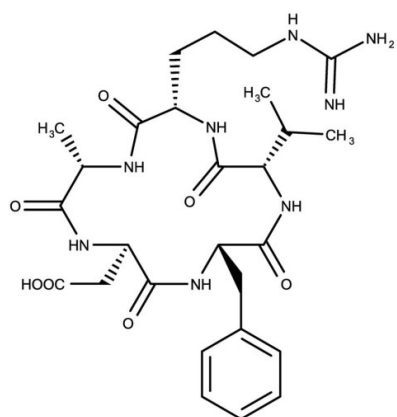
Ordering Information

[Order Online »](#)

BML-AM101-0001	1mg
----------------	-----

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 Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Appearance Solid.

Formula $C_{27}H_{40}N_8O_7$

Formulation Lyophilized.

MW 588.7

Purity ≥98% (HPLC)

Sequence Cyclo[Arg-Ala-Asp-D-Phe-Val]

Solubility Soluble in DMSO (5 mg/ml).

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-
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)