

MMP-3 inhibitor

Inhibitor of matrix metalloproteinase 3 (MMP-3) ($IC_{50}=5\mu M$).

Citations: 2

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-260-165-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

Alternative Name Ac-RCGVPD-NH₂, Stromelysin-1 inhibitor, Matrix metalloproteinase 3 inhibitor

Appearance White to off-white solid.

CAS 158841-76-0

Formula C₂₇H₄₆N₁₀O₉S

MW 686.8

Purity ≥95% (HPLC)

Sequence Ac-Arg-Cys-Gly-Val-Pro-Asp-NH₂

Solubility Soluble in water.

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)