

Z-VRPR-FMK

MALT1 inhibitor

The proteolytic activity of the paracaspase MALT1 is key to NF- κ B and T cell activation. Z-VRPR-FMK is a selective MALT1 inhibitor of such proteolytic activity.

Citations: 10

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-260-166-MC05	0.5mg
ALX-260-166-M001	1mg

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

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

Handling Hygroscopic. Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Z-Val-Arg-Pro-DL-Arg-FMK, MALT1 inhibitor

Appearance White to slightly yellow powder.

Formula $C_{31}H_{49}FN_{10}O_6$

MW 676.8

Purity ≥90% (HPLC)

Sequence Z-Val-Arg-Pro-DL-Arg-fluoromethylketone . TFA

Solubility Soluble in water (1mg/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-
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)