

Furin convertase inhibitor (chloromethylketone)

Highly specific, potent, irreversible, and cell-permeable competitive inhibitor of

proprotein convertases. Highly specific, potent, irreversible, and cell-permeable competitive inhibitor of proprotein convertases: furin/SPC1 ($K_i \sim 1\text{nM}$), SPC2/PC2 ($K_i = 0.36\text{nM}$), SPC3/PC1/PC3 ($K_i = 2.0\text{nM}$), SPC4/PACE4 ($K_i = 3.6\text{nM}$), SPC6/PC5/PC6, and SPC7/LPC/PC7/PC8 ($K_i = 0.12\text{nM}$). Because furin activates viral glycoproteins, this inhibitor is a useful antiviral agent. In addition, it inhibits furin-dependent pro-MMP-2 activation in various cell lines, including U87 and HT1080 cells. Typical concentration range used in tissue culture is 25-100 μM .

Citations: 30

[View Online »](#)

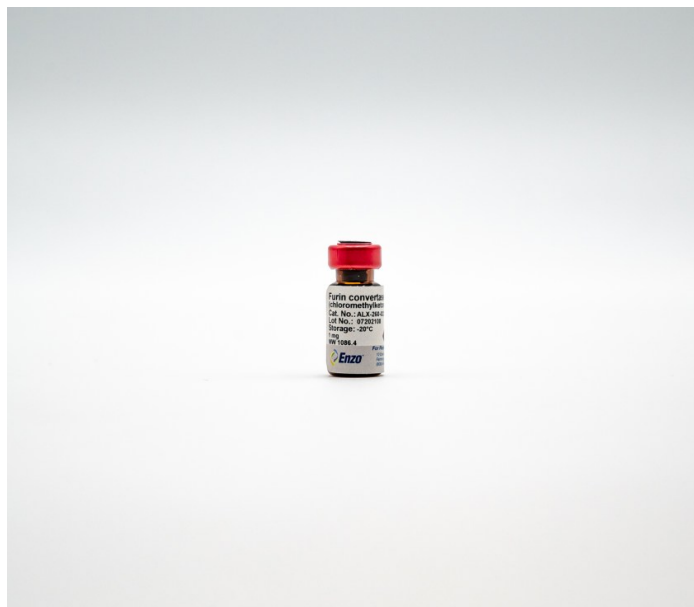
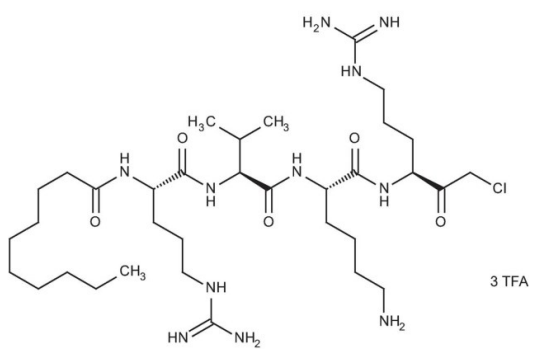
Ordering Information

[Order Online »](#)

ALX-260-022-M001	1mg
ALX-260-022-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 Decanoyl-Arg-Val-Lys-Arg-chloromethylketone, Dec-RVKR-CMK, TACE inhibitor

Appearance White to off-white powder.

Formula $C_{34}H_{66}N_{11}O_5Cl \cdot 3TFA$

MW 744.3 . 342.1

Purity ≥95% (HPLC)

Solubility Soluble in DMSO, methanol or 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)