

Furin convertase substrate (fluorogenic)

Fluorogenic substrate for furin/SPC1/PACE/PCSK3 and other subtilisin/kexin-like proprotein convertases such as Kex2, SPC2/PC2/PCSK2, SPC3/PCSK1, PACE4/SPC4/PCSK6, SPC5/PC4/PCSK4, SPC6/PC5/PC6/PCSK5, and SPC7/LPC/PC7/PC8/PCSK7. Can be used on live cells. Ex: 360-380 nm, Em: 440-460 nm.

Citations: 13

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-260-040-M001	1mg
ALX-260-040-M005	5mg

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Boc-RVRR-AMC, Boc-Arg-Val-Arg-Arg-AMC
Appearance	White to off-white powder.
CAS	136132-77-9
Formula	$C_{38}H_{62}N_{14}O_8$
MW	843.0
Purity	≥95% (HPLC; TLC)
Sequence	Boc-Arg-Val-Arg-Arg-AMC (AMC=7-Amino-4-methylcoumarin)
Solubility	Soluble in water (10mg/ml) or DMSO.

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)