

Boc-Gln-Ala-Arg-AMC

Protease substrate

Fluorogenic substrate for trypsin ($k_{\text{cat}}/K_m = 2.0 \times 10^7 \text{ M}^{-1}\text{sec}^{-1}$; $K_m = 6.0 \mu\text{M}$); matriptase/MT-SP1 and matriptase-2/TMPRSS6; PRSS12; KLK5; KLK11/PRSS20, KLK12, and KLK14; tryptase and thrombin; and dust mite allergen Der p 1. Ex.: 380 nm, Em.: 460 nm, although the following Ex/Em can also be used: 355,375/440,450. This substrate is useful for inhibitor screening and kinetic analysis. Also available: fluorogenic calibration standard AMC (BML-KI144).

Citations: 25

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-P237-0005	5mg
---------------	-----

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Protect from light and moisture.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Boc-QAR-AMC
Appearance	Lyophilized solid.
Formula	$C_{29}H_{42}N_8O_8$
MW	630.7
Purity	≥99% (HPLC)
Sequence	N- <i>tert</i> -butoxycarbonyl-Gln-Ala-Arg-AMC [AMC=7-amino-4-methylcoumarin]
Solubility	Soluble in DMSO (≥10mM) or methanol (50mg/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)