

# Ac-Arg-Leu-Arg-AMC

This fluorogenic peptide is often the substrate of choice for assaying the trypsin-like activity of purified proteasomes. Benefits include a low  $K_m$ , high specific activity, and the fact that the substrate is cleaved exclusively at the R-AMC bond.

Several substrates with R-AMC scissile bonds can be used for assaying trypsin-like activity. One such substrate is Ac-Arg-Leu-Arg-AMC, which benefits from a low  $K_m$ , high specific activity and the fact that it is cleaved exclusively at the R-AMC bond. This fluorogenic peptide is often the substrate of choice for assaying the trypsin-like activity of purified proteasomes.

Citations: 11

[View Online »](#)

## Ordering Information

[Order Online »](#)

BML-AW9785-0005	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** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** Ac-RLR-AMC

**Appearance** Solid.

**Formula**  $C_{30}H_{46}N_{10}O_6$

**Formulation** Lyophilized.

**MW** 642.7

**Purity** ≥97% (HPLC)

**Sequence** Ac-Arg-Leu-Arg-AMC

**Solubility** Soluble in DMSO, methanol or aqueous solutions.



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)