

# Z-LR-AMC

## Cathepsin substrate

Sensitive fluorogenic substrate for cathepsin K ( $K_m=8\mu\text{M}$ ,  $k_{\text{cat}}/K_m=4 \times 10^5 \text{ M}^{-1}\text{s}^{-1}$ ). This substrate is also cleaved by cathepsin B, F, L, L2/V, S. Z-LR-AMC is also cleaved by malaria parasite cysteine proteases falcipain-1, -2, -3, berghepain, and vivapain-2, -3. Ex.: 365nm, Em.: 440nm, although the following Ex/Em can also be used: 355,380/450,460. This substrate is useful for inhibitor screening and kinetic analysis. Also available: fluorogenic calibration standard AMC (BML-KI144).

Citations: 15

[View Online »](#)

## Ordering Information

[Order Online »](#)

BML-P229-0010	10mg
---------------	------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Store, as supplied, at -20°C for up to 1 year. Store solutions at -20°C for up to 3 months.
Handling	Protect from light and moisture.
Long Term Storage	-20°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	Z-Leu-Arg-AMC, Cathepsin K substrate (fluorogenic)
Appearance	White to off-white solid.
Formulation	Lyophilized.
MW	578.7
Purity	≥97% (HPLC)
Sequence	Z-Leu-Arg-AMC . HCl [Z=Cbz=Benzyloxycarbonyl; AMC=7-amino-4-methylcoumarin]
Solubility	Soluble in DMSO to at least 20mM.
Source	Synthetic.

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)