

Suc-Leu-Leu-Val-Tyr-AMC

Proteasome substrate

Fluorogenic substrate for 20S proteasome (Prod. No. BML-PW8720) and other chymotrypsin-like proteases, as well as calpains. Ex.: 380nm Em.: 460nm.

Citations: 105

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-P802-0005	5mg
---------------	-----

Manuals, SDS & CofA

[View Online »](#)

- Fluorogenic substrate of calpains and 20S proteasome
- Used to study the activity of calpains and proteosomal degradation

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	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Suc-LLVY-AMC, Proteasome substrate (fluorogenic), Calpain substrate (fluorogenic)
CAS	94367-21-2
Formula	$C_{40}H_{53}N_5O_{10}$
Formulation	Lyophilized.
Identity	Determined by ESI-MS.
MW	763.9
Purity	≥97% (HPLC)
Sequence	<i>N</i> -Succinyl-Leu-Leu-Val-Tyr-AMC (AMC=7-amino-4-methylcoumarin)
Solubility	Soluble in DMSO (at least 5mM).



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)