

# Ac-IEPD-AMC

## Caspase substrate

Fluorogenic substrate for caspase-8 and granzyme B. Ex.: 360 nm, Em.: 450 nm. The  $K_m$  of recombinant human granzyme B for Ac-IEPD-AMC is 160  $\mu$ M, with a  $V_{max}$  about eight-fold lower than that of Ac-IEPD-pNA (BML-P133).

Citations: 2

[View Online »](#)

## Ordering Information

[Order Online »](#)

|               |     |
|---------------|-----|
| BML-P134-0005 | 5mg |
|---------------|-----|

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

|                   |                                                                                                                                         |
|-------------------|-----------------------------------------------------------------------------------------------------------------------------------------|
| Use/Stability     | As indicated on product label or CoA when stored as recommended. Stock solution in DMSO (e.g. 0.1M) is highly stable at -20°C or -70°C. |
| Handling          | Keep cool and dry.                                                                                                                      |
| Long Term Storage | -20°C                                                                                                                                   |
| Shipping          | Blue Ice                                                                                                                                |

## Regulatory Status

RUO - Research Use Only

## Product Details

|                  |                                                                       |
|------------------|-----------------------------------------------------------------------|
| Alternative Name | Caspase-8 substrate (fluorogenic), Granzyme B substrate (fluorogenic) |
| Appearance       | White to off-white powder.                                            |
| Formula          | $C_{32}H_{41}N_5O_{11}$                                               |
| MW               | 671.8                                                                 |
| Purity           | ≥95%                                                                  |
| Sequence         | N-acetyl-Ile-Glu-Pro-Asp-AMC (7-amino-4-methylcoumarin)               |
| Solubility       | Soluble in 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-  
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)