

# Ac-LEHD-AFC

## Caspase-9 substrate

Fluorogenic substrate for caspase-9.

Citations: 12

[View Online »](#)

## Ordering Information

[Order Online »](#)

|                  |      |
|------------------|------|
| ALX-260-116-M005 | 5mg  |
| ALX-260-116-M010 | 10mg |

Manuals, SDS & CofA

[View Online »](#)

## 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          | Ambient Temperature                                              |

## Regulatory Status

RUO - Research Use Only

## Product Details

|                                |                                                                          |
|--------------------------------|--------------------------------------------------------------------------|
| Alternative Name               | Caspase-9 substrate (fluorogenic)                                        |
| Appearance                     | White to off-white powder.                                               |
| CAS                            | 210345-03-2                                                              |
| Formula                        | $C_{33}H_{38}F_3N_7O_{11}$                                               |
| Formulation                    | Lyophilized.                                                             |
| MW                             | 765.7                                                                    |
| Purity                         | ≥96% (HPLC)                                                              |
| Sequence                       | Ac-Leu-Glu-His-Asp-AFC (AFC=7-Amino-4-trifluoromethylcoumarin)           |
| Solubility                     | Soluble in DMSO.                                                         |
| Technical Info / Product Notes | AFC has an excitation maximum of 400nm and an emission maximum of 505nm. |

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)