

# Ac-LEVD-CHO

Inhibitor of caspase-4 and -5

Reversible inhibitor of caspase-4. Also inhibits caspase-5.

Citations: 3

[View Online »](#)

## Ordering Information

[Order Online »](#)

|                  |     |
|------------------|-----|
| ALX-260-065-M001 | 1mg |
|------------------|-----|

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

|                   |                                                                                                                                       |
|-------------------|---------------------------------------------------------------------------------------------------------------------------------------|
| Use/Stability     | As indicated on product label or CoA when stored as recommended. Stock solutions should be divided into aliquots and stored at -20°C. |
| Handling          | Keep cool and dry.                                                                                                                    |
| Long Term Storage | -20°C                                                                                                                                 |
| Shipping          | Ambient Temperature                                                                                                                   |

## Regulatory Status

RUO - Research Use Only

## Product Details

|                  |                                |
|------------------|--------------------------------|
| Alternative Name | Caspase-4 inhibitor (aldehyde) |
| Appearance       | White powder.                  |
| Formula          | $C_{22}H_{36}N_4O_9$           |
| Formulation      | Lyophilized.                   |
| MW               | 500.6                          |
| Peptide Content  | 70-90%.                        |
| Purity           | ≥95% (HPLC)                    |
| Sequence         | Ac-Leu-Glu-Val-Asp-CHO         |
| Solubility       | Soluble in distilled water.    |

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-  
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)