

# Ac-LEHD-CHO

## Caspase-9 inhibitor

Reversible aldehyde inhibitor of caspase-9 and related cysteine proteases.  
Sequence is from combinatorial experiments.

Citations: 4

[View Online »](#)

### Ordering Information

[Order Online »](#)

|                  |     |
|------------------|-----|
| ALX-260-079-M005 | 5mg |
|------------------|-----|

### Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

|                   |  |
|-------------------|--|
| Use/Stability     | As indicated on product label or CoA when stored as recommended. Keep sealed after removing from the freezer until its temperature equilibrates with room temperature. Dissolve only amount needed at a time to avoid decomposition in solution. |
| Handling          | Keep cool and dry.   |
| Long Term Storage | -20°C  |
| Shipping          | Ambient Temperature  |

## Regulatory Status

RUO - Research Use Only

## Product Details

|                  |   |
|------------------|---|
| Alternative Name | Caspase-9 inhibitor (aldehyde)                                |
| Appearance       | Lyophilized.  |
| Formula          | $C_{23}H_{34}N_6O_9$  |
| MW               | 538.6   |
| Peptide Content  | ≥65%  |
| Purity           | ≥95%  |
| Sequence         | Ac-Leu-Glu-His-Asp-CHO  |
| Solubility       | Soluble in water or DMSO. Further dilute with buffer, pH 7.5. |

Last modified: June 4, 2025



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)