

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. |



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)