

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 »](#)

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

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com