

Ac-LEHD-AMC

Caspase-9 substrate

Fluorogenic substrate for caspase-9 and related cysteine proteases. The LEHD sequence is derived from combinatorial experiments. Ex.:340-360nm, Em.:440-460nm.

Citations: 13

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-260-080-M001	1mg
ALX-260-080-M005	5mg

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
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 substrate (fluorogenic)
Appearance	White to off-white powder.
CAS	292633-16-0
Formula	$C_{33}H_{41}N_7O_{11}$
MW	711.7
Peptide Content	75-95%
Purity	≥98% (HPLC)
Sequence	Ac-Leu-Glu-His-Asp-AMC (AMC=7-Amino-4-methylcoumarin)
Solubility	Soluble in water, DMSO or phosphate buffer, pH 7.5.
Technical Info / Product Notes	AMC has an excitation maximum of 340-360nm and an emission maximum of 440-460nm.

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)