

Ac-LEVD-AFC

Caspase-4 substrate

Fluorogenic substrate for caspase-4. Similar to the fluorogenic substrate Ac-LEVD-AMC (Prod. No. ALX-260-083); the AFC fluorophore is more sensitive.

Ordering Information

[Order Online »](#)

ALX-260-084-M001	1mg
ALX-260-084-M005	5mg

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Dissolve only as much as needed at one time because of the possibility of decomposition in solution. Store solution at -20°C. Protect from condensation: when taking the sample out of the container, wait until it warms up to room temperature, then open it.
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 substrate (fluorogenic)
Appearance	White to off-white powder.
Formula	$C_{32}H_{40}F_3N_5O_{11}$
MW	727.7
Peptide Content	75-95%.
Purity	≥95% (HPLC)
Sequence	Ac-Leu-Glu-Val-Asp-AFC (AFC=7-Amido-4-trifluoromethylcoumarin)
Solubility	Soluble in DMSO; dilute with distilled water or buffer; also soluble in phosphate buffer, pH 7.5-8.0.
Technical Info / Product Notes	AFC has an excitation maximum of 400nm and an emission maximum of 505nm.

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-
de@enzolifesciences.com](mailto:info-de@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)