

Ac-WEHD-AFC

Substrate for caspase-1, -4-, -5 and -14

Fluorogenic substrate for caspase-1, -4, -5 and -14.

Citations: 3

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-260-117-M005	5mg
ALX-260-117-M010	10mg

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Solution in DMSO is stable for 6-8 months when stored at -20°C.
Handling	Keep cool and dry.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Caspase-1 substrate (fluorogenic), Caspase-5 substrate (fluorogenic)
Appearance	White to yellow powder.
Formula	$C_{38}H_{37}F_3N_8O_{11}$
MW	838.8
Peptide Content	70-90%
Purity	≥98% (HPLC)
Sequence	Ac-Trp-Glu-His-Asp-AFC (AFC=7-Amino-4-trifluoromethylcoumarin)
Solubility	Soluble in DMSO.
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-
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)