

Ac-WEHD-AMC

Substrate for caspase-1, -4 and -5

Fluorogenic substrate for caspase-1 ($K_m=4\mu\text{M}$), -4 ($K_m=31\mu\text{M}$);and -5 ($K_m=15\mu\text{M}$). This fluorogenic caspase-1 substrate with an optimal tetrapeptide sequence is cleaved 50-fold more efficiently than Ac-YVAD-AMC (Prod. No. ALX-260-024).

Citations: 7

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-260-057-M001	1mg
ALX-260-057-M005	5mg

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Handling Protect from light and moisture.

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), Caspase-4 substrate (fluorogenic)

Formula $C_{38}H_{40}N_8O_{11}$

Formulation Lyophilized.

MW 784.8

Peptide Content 70-90%

Purity ≥98% (HPLC)

Sequence Ac-Trp-Glu-His-Asp-AMC (AMC=7-Amino-4-methylcoumarin)

Solubility Soluble in DMSO (100mM, 75.5mg/ml). Dilute with distilled water or buffer.

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)