

Ac-VEID-AMC

Caspase-6 substrate

Fluorogenic substrate for caspase-6 ($K_m=30\mu\text{M}$, $K_{cat}/K_m=168000\text{M}^{-1}\text{se}^{-1}$) and related cysteine proteases. Sequence is based on lamin A site of cleavage. Ex.:340-360nm, Em.:440-460nm.

Citations: 10

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-260-064-M001	1mg
ALX-260-064-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. Keep cool and dry.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Caspase-6 substrate (fluorogenic)

Appearance White to off-white powder.

CAS 219137-97-0

Formula $C_{32}H_{43}N_5O_{11}$

MW 673.7

Peptide Content 70-90%

Purity ≥95% (HPLC)

Sequence Ac-Val-Glu-Ile-Asp-AMC (AMC=7-Amino-4-methylcoumarin)

Solubility Soluble in DMSO; dilute with distilled water or buffer; also soluble in 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-
be@enzolifesciences.com](mailto:info-be@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)