

Ac-VEID-AFC

Caspase-6 inhibitor

Fluorogenic substrate for caspase-6 ($K_m=30\mu\text{M}$; $K_{cat}/K_m=168000\text{ M}^{-1}\text{sec}^{-1}$) and related cysteine proteases. Sequence is based on lamin A site of cleavage.

Citations: 3

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-260-111-M005	5mg
ALX-260-111-M010	10mg

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Caspase-6 substrate (fluorogenic)
Formula	$C_{32}H_{40}F_3N_5O_{11}$
MW	727.7
Purity	≥98% (HPLC)
Sequence	Ac-Val-Glu-Ile-Asp-AFC (AFC=7-Amino-4-trifluoromethylcoumarin)
Solubility	Soluble in DMSO or dimethyl formamide.
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

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com