

Ac-YVAD-AFC

Substrate for caspase-1 and -4

Fluorogenic substrate for caspase-1 and -4.

Citations: 3

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-260-108-M005	5mg
ALX-260-108-M010	10mg

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Solid product is stable for 1 year in desiccator when stored at room temperature. Stock solutions are stable for 6 – 8 months when stored at –20°C. Keep sealed after removing from the freezer until its temperature equilibrates with room temperature.
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-4 substrate (fluorogenic)
Appearance	White to off-white powder.
Formula	$C_{33}H_{36}F_3N_5O_{10}$
MW	719.7
Purity	≥97% (HPLC)
Sequence	Ac-Tyr-Val-Ala-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.



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)