

# Z-YVAD-AFC

## Caspase-1 substrate

Fluorogenic substrate for caspase-1. AFC has an excitation maximum of 400nm and an emission maximum of 505nm. This product is similar to Ac-YVAD-AMC, Prod. No. ALX-260-024, but the AFC fluorophore has a greater Stokes' shift upon cleavage and the N-terminal is blocked by a benzyloxycarbonyl moiety, which increases cell permeability.

Citations: 12

[View Online »](#)

## Ordering Information

[Order Online »](#)

|                  |     |
|------------------|-----|
| ALX-260-035-M001 | 1mg |
| ALX-260-035-M005 | 5mg |

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

|                   |                                                                                                                     |
|-------------------|---------------------------------------------------------------------------------------------------------------------|
| Use/Stability     | As indicated on product label or CoA when stored as recommended. DMF or DMSO stock solution highly stable at -20°C. |
| Handling          | Protect from light.                                                                                                 |
| Long Term Storage | -20°C                                                                                                               |
| Shipping          | Ambient Temperature                                                                                                 |

## Regulatory Status

RUO - Research Use Only

## Product Details

|                  |                                                                 |
|------------------|-----------------------------------------------------------------|
| Alternative Name | Caspase-1 substrate (fluorogenic)                               |
| Appearance       | White to off-white powder.                                      |
| CAS              | 201608-13-1                                                     |
| Formula          | $C_{39}H_{40}N_5O_{11}F_3$                                      |
| MW               | 811.8                                                           |
| Purity           | ≥95% (HPLC).Single spot (TLC)                                   |
| Sequence         | Z-Tyr-Val-Ala-Asp-AFC (AFC = 7-Amido-4-trifluoromethylcoumarin) |
| Solubility       | Soluble in DMSO, DMF or methanol; insoluble in water.           |

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)