

Ac-DEVD-AFC

Caspase-3 substrate

Fluorogenic substrate for caspase-3 (CPP32), with a $K_m=9.7\mu M$ and related cysteine proteases. Sequence is based on PARP cleavage at Asp²¹⁶ for caspase-3. Similar to Ac-DEVD-AMC (Prod. No. ALX-260-031) but the AFC fluorophore has a greater Stokes' shift upon cleavage than AMC. Reaction can be monitored quantitatively or visually using a hand-held long-UV lamp and visualizing a blue to green shift in fluorescence upon cleavage. Ex.: 400nm, Em.: 505nm.

Citations: 32

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-260-032-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-3 substrate (fluorogenic)
Appearance	White to off-white powder.
CAS	201608-14-2
Formula	$C_{30}H_{34}F_3N_5O_{13}$
MW	729.6
Peptide Content	65-95%
Purity	≥96% (HPLC)
Sequence	Ac-Asp-Glu-Val-Asp-AFC (AFC = 7-Amino-4-trifluoromethylcoumarin)
Solubility	Soluble in dimethyl formamide, DMSO or methanol; slightly soluble in water (0.4mg/ml).
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](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-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)

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