

Ac-VDVAD-AFC

Caspase-2 substrate

Fluorogenic substrate for caspase-2; similar to Ac-VDVAD-AMC but the AFC fluorophore has a greater Stokes' shift upon cleavage. VDVAD has been found to be a preferred cleavage site for caspase-2 (Ich-1L). Increases in Ac-VDVAD-AFC cleavage correlated with losses in procaspase-2 (conversion to active caspase-2) in human neuroblastoma lines induced to apoptosis with C2-ceramide and NO.

Ordering Information

[Order Online »](#)

ALX-260-112-M005	5mg
ALX-260-112-M010	10mg

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Warm up to room temperature before opening. Keep cool and dry.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	Caspase-2 substrate (fluorogenic)
Appearance	White to off-white powder.
Formula	$C_{33}H_{41}F_3N_6O_{12}$
MW	770.7
Peptide Content	75-95%
Purity	≥98% (HPLC)
Sequence	Ac-Val-Asp-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

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-fr@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