

Ac-IETD-AMC

Caspase and granzyme substrate

Fluorogenic substrate for caspase-6 and -8 and granzyme B. Also a substrate for caspase-10.

Citations: 12

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-260-042-M001	1mg
------------------	-----

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	Granzyme B substrate (fluorogenic), Caspase-8 substrate (fluorogenic)
Appearance	White to off-white powder.
Formula	$C_{31}H_{41}N_5O_{12}$
MW	675.7
Peptide Content	70-90%
Purity	≥97% (HPLC)
Sequence	Ac-Ile-Glu-Thr-Asp-AMC (AMC=7-Amino-4-methylcoumarin)
Solubility	Soluble in DMSO. Dilute with buffer, pH 7.5.
Technical Info / Product Notes	AMC has an excitation maximum of 340-360nm and an emission maximum of 440-460nm.

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-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