

Ac-IETD-AFC

Caspase and granzyme substrate

Fluorogenic substrate for caspase-8/caspase-3 processing enzyme and granzyme B. Also a substrate for caspase-10. The IETD sequence is based on caspase-3 proenzyme cleavage site.

Citations: 11

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-260-110-M005	5mg
ALX-260-110-M010	10mg

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Handling Keep cool and dry.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Caspase-8 substrate (fluorogenic), Granzyme B substrate (fluorogenic)

Appearance White to off-white powder.

Formula $C_{31}H_{38}F_3N_5O_{12}$

MW 729.7

Purity ≥98% (HPLC)

Sequence Ac-Ile-Glu-Thr-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.

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