

Ac-LETD-AFC

Caspase-8 substrate

Fluorogenic substrate for caspase-8.

Citations: 4

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. 20mM stock solution in DMSO is stable at -20°C for 6-8 months. When taking out of the freezer wait until it warms up to room temperature, then open it.
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)
Appearance	White to off-white powder.
CAS	210345-02-1
Formula	$C_{31}H_{38}F_3N_5O_{12}$
MW	729.7
Purity	≥97% (HPLC)
Sequence	Ac-Leu-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](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)