

Ac-IETD-CHO

Caspase and granzyme inhibitor

Reversible inhibitor of caspase-6, -8 and granzyme B. Also inhibits caspase-10.

Citations: 3

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-260-043-M001	1mg
ALX-260-043-M005	5mg

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Use always fresh solutions.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Caspase-8 inhibitor (aldehyde), Granzyme B inhibitor (aldehyde)
CAS	191338-86-0
Formula	$C_{21}H_{34}N_4O_{10}$
Formulation	Lyophilized.
MW	502.5
Peptide Content	70-90%
Purity	≥96% (HPLC)
Sequence	Ac-Ile-Glu-Thr-Asp-CHO
Solubility	Soluble in distilled water at 10 mg/mL and in DMSO at 1 mg/mL.

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-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

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

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