

Caspase fluorometric substrate set

Fluorometric substrates for assaying activities of members of caspase family proteases. All substrates are provided in liquid *ready-to-use* form.

Ordering Information

[Order Online »](#)

ALX-850-229-KI01	1Set
------------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application	Fluorescence microscopy
Contents	Contains 125µl (1mM) each of the following substrates: – Caspase-1 Substrate, Ac-YVAD-AFC – Caspase-2 Substrate, Ac-VDVAD-AFC – Caspase-3 Substrate, Ac-DEVD-AFC – Caspase-5 Substrate, Ac-WEHD-AFC – Caspase-6 Substrate, Ac-VEID-AFC – Caspase-8 Substrate, Ac-IETD-AFC – Caspase-9 Substrate, Ac-LEHD-AFC
Formulation	Liquid. <i>Ready-to-use</i> in DMSO.

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-
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)