

Caspase-2

fluorometric assay kit

Based on the detection of AFC (7-amino-4-trifluoromethyl coumarin) after cleavage from the labelled substrate VDVAD-AFC.

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-850-214-KI01	25 tests
------------------	----------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	Store Cell Lysis Buffer and 2X Reaction Buffer at 4°C after opening.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application	Activity assay, Fluorescent detection
Contents	25ml Cell Lysis Buffer, 2ml 2X Reaction Buffer, 125µl VDVAD-AFC Substrate (1mM), 100µl DTT.
Technical Info / Product Notes	VDVAD-AFC emits blue light ($\lambda_{\text{max}} = 400\text{nm}$); upon cleavage of the substrate by caspase-2 or related caspases, free AFC emits a yellow-green fluorescence ($\lambda_{\text{max}} = 505\text{nm}$), which can be quantified using a fluorometer or a fluorescence microtiter plate reader. Comparison of the fluorescence of AFC from an apoptotic sample with an uninduced control allows determination of the fold increase in caspase-2 activity. Biocompare Article: Keep an Eye on Apoptosis with Caspase Assays
UniProt ID	P42575

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-be@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