

# Caspase-3 fluorometric assay kit

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

Citations: 8

[View Online »](#)

## Ordering Information

[Order Online »](#)

ALX-850-216-KI01	25 tests
------------------	----------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

**Handling** Store Lysis Buffer and 2X Reaction Buffer at 4°C after opening. Protect from light.

**Long Term Storage** -20°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** CPP32, Yama, Apopain

**Application** Activity assay, Fluorescent detection

**Contents** 25ml Cell Lysis Buffer, 2ml 2X Reaction Buffer, 125µl DEVD-AFC (1mM), 100µl DTT (1M).

**Technical Info / Product Notes** DEVD-AFC emits blue light ( $\lambda_{\text{max}} = 400\text{nm}$ ); upon cleavage of the substrate by caspase-3 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-3 activity.

Biocompare Article: [Keep an Eye on Apoptosis with Caspase Assays](#)

**UniProt ID** P42574

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