

Caspase-1 assay kit for drug discovery

NOW INCLUDES BOTH COLORIMETRIC AND FLUOROGENIC SUBSTRATES! This kit is a complete assay system to measure protease activity of recombinant caspase-1. Cleavage of a tetrapeptide substrate is monitored colorimetrically at 405 nm or fluorimetrically (Ex. 360 nm/Em. 460 nm). Assays are performed in a convenient 96-well plate format.

Citations: 2

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-AK701-0001	96 wells
----------------	----------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application

Activity assay, Colorimetric detection, Fluorescent detection, HTS

Compatibility

This product is compatible with the [Absorbance 96 Plate Reader](#).

Contents

Caspase-1 enzyme (human) (recombinant), Ac-YVAD-pNA colorimetric substrate, Ac-YVAD-AMC fluorogenic substrate, calibration standards, aldehyde inhibitor, assay buffer, 1/2-volume 96-microwell plate, and instruction booklet.

Technical Info / Product Notes

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

UniProt ID

P29466

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)