

Caspase-3 cell assay kit

NOW INCLUDES BOTH COLORIMETRIC AND FLUOROGENIC SUBSTRATES! This kit is a complete assay system for measuring caspase-3 and caspase-3-like activity (DEVDase) in cellular extracts. It is performed in a convenient 96-well format with all reagents necessary for preparing cell extracts, measuring caspase activity and calibrating the assay. *PLUS*, Caspase-3 enzyme is included! The caspase-3, is useful as a positive control, for testing cellular extracts for endogenous inhibitors, and more.

Citations: 10

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-AK703-0001	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

Alternative Name CPP32, Yama, Apopain

Application Activity assay, Colorimetric detection, Fluorescent detection, HTS

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

Contents

Caspase-3 enzyme (human, recombinant) (Prod. No. BML-SE169-9090) (300U, (30µl); 10 u/µl in assay buffer. One U=1 pmol/min at 30°C) Storage: -70°C, avoid freeze/thaw cycles **pNA Substrate** (Ac-DEVD-pNA, MW=637) (Prod. No. BML-P412-9090) (1ml; 2mM (1.3 mg/ml) in assay buffer) Storage: -70°C **pNA calibration standard** (p-nitroaniline; MW=138) (Prod. No. BML-KI106) (1 ml, 50µM in assay buffer. A405nm=0.525) Storage: -70°C **AMC Substrate** (Ac-DEVD-AMC; MW=676) (Prod. No. BML-P411-9090) (1 ml; 0.3mM (0.20 mg/ml) in assay buffer Storage: -70°C **AMC calibration standard** (7-amino-4-methylcoumarin; MW=175)(Prod. No. BML-KI107) (1 ml; 30µM in assay buffer) Storage: -70°C **Inhibitor** (Ac-DEVD-CHO; MW=502) (Prod. No. BML-P410-9090) (50µl; 0.1mM (0.05 mg/ml) in DMSO) Storage: -70°C **Assay Buffer** (Prod. No. BML-KI-111) (20 ml; 50mM HEPES, pH 7.4, 100mM sodium chloride, 0.1% CHAPS, 10mM dithiothreitol, 1mM ethylenediaminetetraacetic acid, 10% glycerol) Storage: -70°C **Cell lysis buffer** (50mM HEPES, pH 7.4, 0.1% CHAPS, 5mM dithiothreitol, 0.1mM ethylenediaminetetraacetic acid) (Prod. No. BML-KI117) (30 ml; liquid in white screw-cap plastic bottle) Storage: -70°C **½-Volume microplate** (Prod. No. BML-KI101) Storage: Room temperature

Technical Info / Product Notes

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

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)