

CV-Caspase 3 & 7 detection kit

A consistent feature of mammalian cell apoptosis is the activation of caspases 3 and/or 7. Based on the dye cresyl violet and the caspase 3/7 DEVD peptide recognition sequence, this kit allows the detection of caspases 3 and 7 in living cells by fluorescence microscopy. When cleaved by caspase 3 or 7, the mono and unsubstituted cresyl violet fluoresce when excited at 550-590 nm. Once cells are grown, and apoptosis induced, staining is completed in just a few minutes. Hoechst 33342 and Acridine Orange are also included to allow visualization of nuclei and lysosomes.

Citations: 4

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-AK118-0001	1Kit
----------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling	Store the unopened kit and each unopened component at 2-8°C until the expiration date. Once reconstituted with DMSO, use CR(DEVD) ₂ immediately, or store at ≤-20°C up to 6 months, protected from light and thawed no more than twice during that time.
Long Term Storage	+4°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Application	Fluorescence microscopy, Fluorescent detection
Contents	CR(DEVD) ₂ reagent, Acridine orange, Hoechst 33342 stain.

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)