

CV-Cathepsin L detection kit

Elevated cathepsin enzyme activity often indicates the presence of a pathological condition including cancer, Alzheimer's disease, autoimmune disease, and osteoporosis. Based on the dye cresyl violet and the cathepsin L PheArg peptide recognition sequence, this kit allows the detection of cathepsin L activity in living cells by fluorescence microscopy. Once cells are grown, staining is completed in just a few minutes. Hoechst 33342 stain and Acridine Orange are also included to allow visualization of nuclei and lysosomes.

Citations: 1

[View Online »](#)

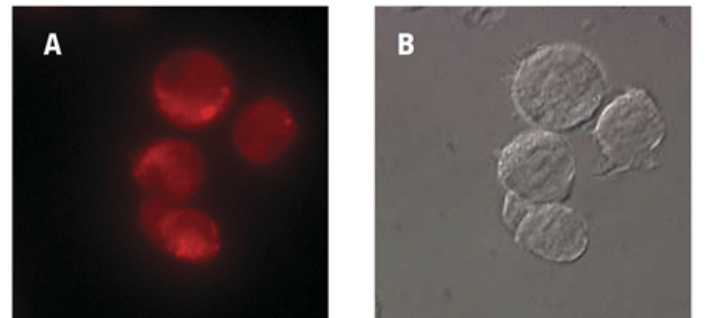
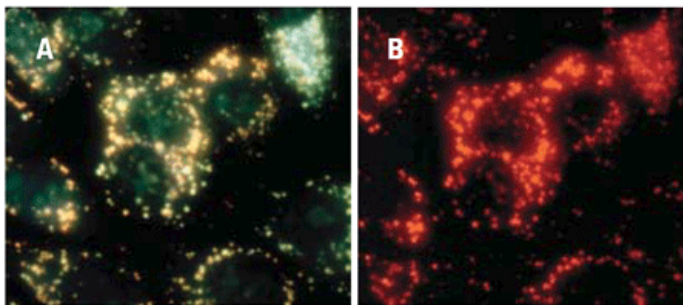
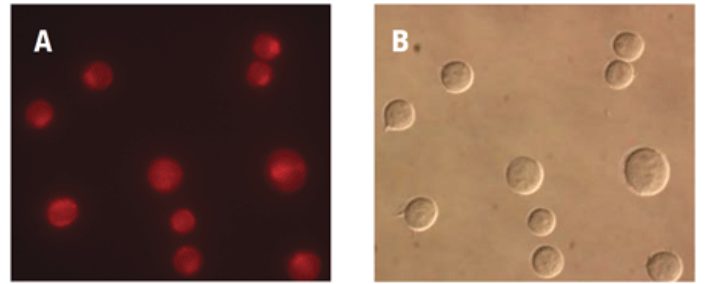
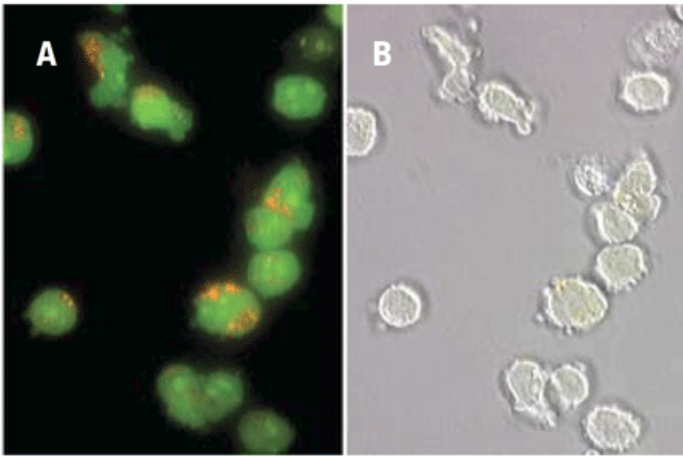
Ordering Information

[Order Online »](#)

BML-AK127-0001	1Kit
----------------	------

Manuals, SDS & CofA

[View Online »](#)



Long Term Storage +4°C

Regulatory Status RUO - Research Use Only

Application	Fluorescence microscopy, Fluorescent detection
--------------------	--

Contents CV-(FR)₂ reagent, Acridine Orange, Hoechst 33342 stain.

UniProt ID P07711 (cathepsin L1), O60911 (cathepsin L2)



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@enzolifesciences.com)