

# CV-Cathepsin B 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 B ArgArg peptide recognition sequence, this kit allows the detection of cathepsin B 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: 10

[View Online »](#)

## Ordering Information

[Order Online »](#)

BML-AK125-0001	1Kit
----------------	------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Application** Fluorescence microscopy, Fluorescent detection

**Application Notes** Cited sample types include mouse microglial cells, human monocytic leukemia cells, and human PBMC.

**Contents** CV-(RR)<sub>2</sub> reagent, Acridine Orange, Hoechst 33342 stain

**UniProt ID** P07858



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)