## 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** 

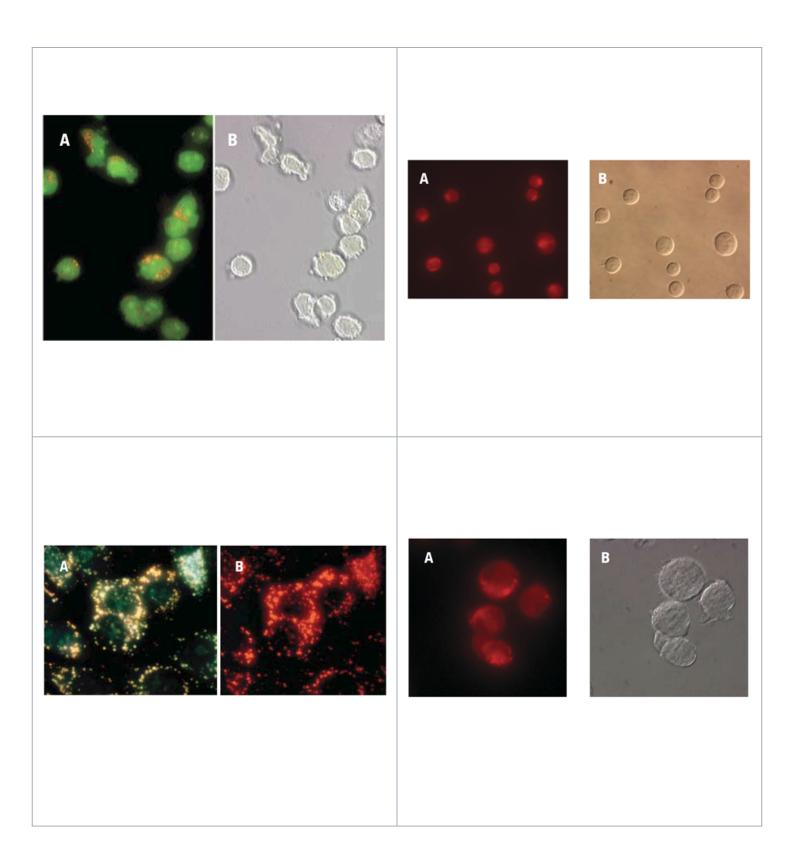
**Order Online** »

BML-AK127-0001

1Kit

Manuals, SDS & CofA

View Online »



## **Handling & Storage**

**Long Term Storage** +4°C

**Shipping** Blue Ice

Regulatory Status RUO - Research Use Only

## **Product Details**

**Application** Fluorescence microscopy, Fluorescent detection

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

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

Last modified: May 29, 2024



info-