CV-Cathepsin K 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 K LeuArg peptide recognition sequence, this kit allows the detection of cathepsin K 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.

Ordering Information

Order Online »

BML-AK126-0001

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-(LR)₂ reagent, Acridine Orange, Hoechst 33342 stain.

P43235 **UniProt ID**

Last modified: May 29, 2024



info-

eu@enzolifesciences.com