

# Cathepsin D-E Activity Assay

**A plate based fluorometric assay to measure Cathepsin D activity in a variety of biological samples with a sensitivity able to detect Cathepsin B in 500 cells or less.**

Cathepsin D performs a variety of functions in the cell, including recycling of protein through autophagy, maturation of proteins by cleaving precursor proteins and apoptosis. Cathepsin D has a dual role with respect to apoptosis. On the one hand, under conditions of cellular stress such as DNA damage or oxidative stress, Cathepsin D is released from the lysosomes and activates apoptotic pathways by cleaving pro-apoptotic proteins. On the other hand, Cathepsin D can prevent apoptosis by regulating the stability and activity of anti-apoptotic proteins, inhibiting apoptotic pathways, enhancing cellular survival and tissue protection. AkrisBio's Cathepsin D Activity Assay is a simple, sensitive means of measuring Cathepsin D activity in a variety of biological samples with a sensitivity able to detect Cathepsin B in 500 cells or less.

## Ordering Information

[Order Online »](#)

AKR-MA-0154	100 wells
-------------	-----------

## Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage      -20°C

Shipping                      Blue Ice

**Regulatory Status**      RUO - Research Use Only

## Product Details

Application                      Fluorescent detection

Species Reactivity              Species independent

**Technical Info / Product Notes**      For the Original Manufacturer's data sheet please [click here](#).



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)