

Cathepsin D Inhibitor Assay Kit ab126779

[1 References](#) [1 Image](#)

Overview

| | |
|-------------------------|--|
| Product name | Cathepsin D Inhibitor Assay Kit |
| Detection method | Fluorescent |
| Sample type | Inhibitor compounds |
| Assay type | Quantitative |
| Assay time | 1h 00m |
| Product overview | Cathepsin D Inhibitor Screening Kit (ab126779) is a fluorescence-based assay that utilizes the preferred cathepsin D substrate sequence GKPILFFRLK(Dnp)-D-R-NH ₂) labeled with MCA. Cathepsin D will cleave the synthetic substrate to release the quenched fluorescent group MCA, which can then easily be measured using a fluorometer or fluorescence plate reader at Ex/Em = 328/460 nm. The relative efficacy of test inhibitors are compared to the positive control inhibitor, Pepstatin A (IC ₅₀ < 0.1 nM). The Cathepsin D assay is simple, straightforward, and can be adapted to 96-well plate assays and is suitable for high throughput screening (HTS). Visit our FAQs page for tips and troubleshooting. |

| | |
|--------------|---|
| Notes | <p>This product is manufactured by BioVision, an Abcam company and was previously called K148 Cathepsin D Inhibitor Screening Kit (Fluorometric). K148-100 is the same size as the 100 test size of ab126779.</p> <p>Apoptosis can be mediated by mechanisms other than the traditional caspase-mediated cleavage cascade. There is growing recognition that alternative proteolytic enzymes such as the lysosomal cathepsin proteases may initiate or propagate pro-apoptotic signals. Cathepsins are lysosomal enzymes that are also used as sensitive markers in various toxicological investigations.</p> |
|--------------|---|

| | |
|-----------------|-------------------|
| Platform | Microplate reader |
|-----------------|-------------------|

Properties

| | |
|-----------------------------|----------------------------|
| Storage instructions | Please refer to protocols. |
|-----------------------------|----------------------------|

| Components | 100 tests |
|--------------------|-----------|
| CD Reaction Buffer | 1 x 10ml |
| Human Cathepsin D | 1 vial |

| Components | 100 tests |
|----------------------|-----------|
| Pepstatin A Solution | 1 x 20µl |
| Substrate II | 1 x 0.2ml |

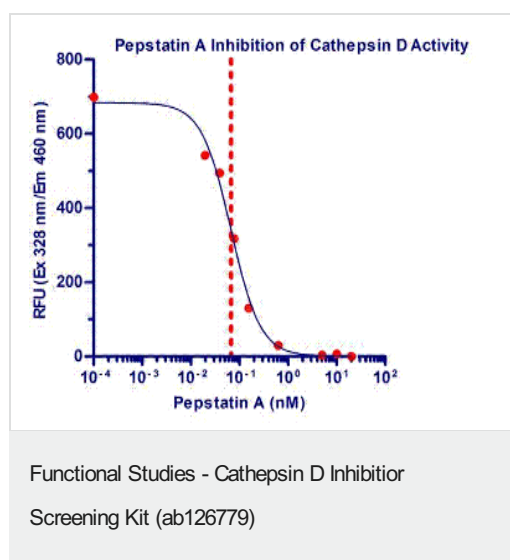
Relevance

Cathepsin D is a normal lysosomal protease that is expressed in all cells. It is an aspartyl protease with a pH optimum in the range of 3-5, and contains two N linked oligosaccharides. Cathepsin D is synthesized as an inactive pro enzyme. Activation involves the proteolytic removal of the 43 amino acid profragment and an internal cleavage to generate the two chain form made up of 34 and 14 kDa subunits. Cathepsin D contains the mannose-6-phosphate lysosomal localization signal that targets the enzyme to the lysosomal compartment where it functions in the normal degradation of proteins. In certain tumor cells, Cathepsin D is abnormally processed and is secreted in its precursor form. Numerous clinical studies as well as in vitro evidence suggest that cathepsin D plays an important role in malignant transformation and may be a useful prognostic indicator for breast cancer and possibly Alzheimer's disease.

Cellular localization

Lysosome. Melanosome. Note=Identified by mass spectrometry in melanosome fractions from stage I to stage IV.

Images



Typical Pepstatin A Inhibition Profile of Cathepsin D Activity. The red line denotes an IC_{50} value of 0.067 nM. Results were analyzed by fluorescence plate reader according to the kit instructions.

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors