

Product datasheet

Cathepsin S Activity Assay Kit (Fluorometric) ab65307

3 References 2 Images

Overview

Product name	Cathepsin S Activity Assay Kit (Fluorometric)
Sample type	Tissue Extracts, Cell Lysate
Assay type	Enzyme activity
Assay time	2h 00m
Product overview	<p>Cathepsin S Activity Assay Kit (ab65307) is a fluorescence-based assay that utilizes the preferred cathepsin-S substrate sequence VVR labeled with AFC (amino-4-trifluoromethyl coumarin). Cell lysates or other samples that contain cathepsin-S will cleave the synthetic substrate VVR-AFC to release free AFC. The released AFC can easily be quantified using a fluorometer or fluorescence plate reader.</p> <p>Visit our FAQs page for tips and troubleshooting.</p>

Notes

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 proapoptotic signals. Cathepsins are lysosomal enzymes that are also used as sensitive markers in various toxicological investigations.

Tested applications **Suitable for:** Functional Studies

Properties

Storage instructions Store at -20°C. Please refer to protocols.

Components	Identifier	100 tests
CS Cell Lysis Buffer	WM	1 x 25ml
CS Inhibitor (1 mM)	Red	1 x 20µl
CS Reaction Buffer	NM	1 x 5ml
CS Substrate Ac-VVR-AFC (10 mM)	Brown	1 x 200µl

Relevance

The cathepsin family of proteases contains several diverse classes of enzymes. The cysteine protease class comprises cathepsin B, H, K, L, O, and S. The aspartyl protease class contains cathepsin D and E. Cathepsin G belongs to the serine protease class. Cathepsins are involved

in various cellular events such as peptide biosynthesis, protein degradation, and apoptosis. Cathepsin S has been shown to be able to function as an elastase over a broad pH range in alveolar macrophages.

Cellular localization

Lysosome

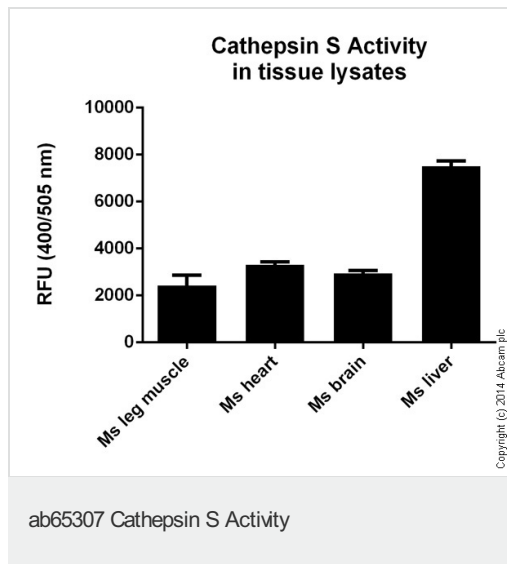
Applications

Our [Abpromise guarantee](#) covers the use of **ab65307** in the following tested applications.

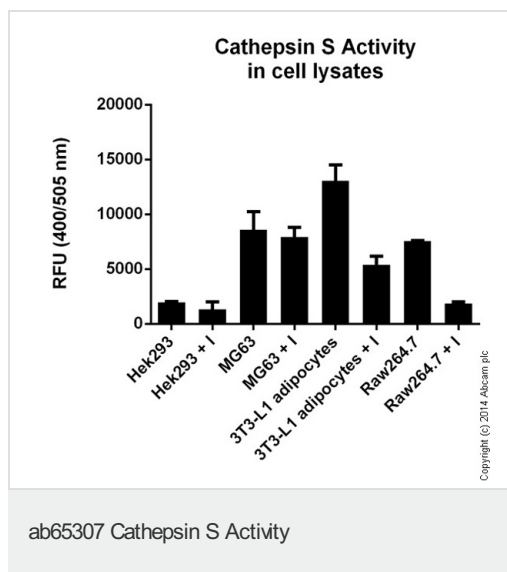
The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

Application	Abreviews	Notes
Functional Studies		Use at an assay dependent dilution.

Images



Cathepsin S Activity measured in tissue lysates (each with protein concentration of 4 mg/mL) after 1 hour of incubation.



Cathepsin S Activity measured in cell lysates (each with cells concentration of 4×10^7 cells/mL) after 1 hour of incubation with and without inhibitor (I).

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 <http://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