

Technical Bulletin

Cathepsin B Activity Assay Kit

Catalog Number MAK387

Product Description

Apoptosis can be mediated by mechanisms other than the traditional caspase-mediated cleavage cascade. There is growing recognition that alternative proteolytic enzymes such as lysosomal cathepsin proteases may initiate or propagate proapoptotic signals. Cathepsins are lysosomal enzymes that are also used as sensitive markers in various toxicological investigations.

The Cathepsin B Activity Assay kit uses a fluorescence-based assay that utilizes the preferred cathepsin B substrate sequence RR labeled with AFC (amino-4-trifluoromethyl coumarin). Cell lysates or other samples that contain cathepsin B cleave the synthetic substrate RR-AFC to release free AFC. The released AFC can easily be quantified fluorometrically. The cathepsin B assay is simple and straightforward. Assay conditions have been optimized to obtain the maximal activity.

The kit is suitable for the detection of Cathepsin B activity in cell lysates.

Components

The kit is sufficient for 100 fluorometric assays in 96-well plates.

- | | |
|---------------------------------------------------------------|--------|
| • CB Cell Lysis Buffer
Catalog Number MAK387A | 25 mL |
| • CB Reaction Buffer
Catalog Number MAK387B | 5 mL |
| • CB Substrate Ac-RR-AFC
(10 mM)
Catalog Number MAK387C | 200 µL |
| • CB Inhibitor (1 mM)
Catalog Number MAK387D | 20 µL |

Reagents and Equipment
Required but Not Provided

- Pipetting devices and accessories (e.g., multichannel pipettor)
- Fluorescence multiwell plate reader
- Flat bottom, opaque, white or black 96-well plate for enhanced sensitivity. Cell culture or tissue culture treated plates are **not** recommended.
- Microcentrifuge
- 7-Amino-4-(trifluoromethyl)coumarin (AMC) (Catalog Number 248924) (optional)

Precautions and Disclaimer

For R&D use only. Not for drug, household, or other uses. Please consult the Safety Data Sheet for information regarding hazards and safe handling practices.

Storage/Stability

The kit is shipped on wet ice. Store components at -20 °C.

Preparation Instructions

Briefly centrifuge small vials prior to opening.

CB Cell Lysis Buffer and CB Reaction Buffer:
Store at 2-8 °C after opening.

CB Substrate Ac-RR-AFC (10 mM): Protect from light.

Procedure

Sample Preparation

1. Induce sample cells to increase Cathepsin B activity by adding test chemical. For unknown test chemicals, it is recommended to prepare several cell cultures and add different concentrations of test chemical.
2. Prepare a separate cell culture with **no** test chemical inducer for use as an Uninduced Control.
3. Collect cells ($1-5 \times 10^6$ cells) by centrifugation.
4. Lyse cells in 50 μ L of chilled CB Cell Lysis Buffer.
5. Incubate cells on ice for 10 minutes.
6. Centrifuge at top speed in a microcentrifuge for 5 minutes.
7. Transfer the supernatants to new tubes.
8. Add 50 μ L of each cell lysate, including Uninduced Control cell lysate, to a 96-well plate. Note: If not running multiple sample cell lysates with different amounts of test chemical, use 50-200 μ g of cell lysates (diluted to a total volume of 50 μ L of CB Cell Lysis Buffer) if protein concentration has been measured.

Negative Cell Lysate Control (Optional)

Add 50 μ L of Uninduced Control cell lysate and 2 μ L of CB Inhibitor (1 mM) to designated well.

Assay Procedure

1. Add 50 μ L of CB Reaction Buffer to all Induced Sample and Control (Uninduced and optional Negative Cell Lysate) wells.
2. Add 2 μ L of the 10 mM CB Substrate Ac-RR-AFC (200 μ M final concentration) to each well.
3. Incubate at 37 °C for 1-2 hours.

Measurement

Measure the fluorescence (RFU) at $\lambda_{Ex} = 400 \text{ nm}/\lambda_{Em} = 505 \text{ nm}$.

Results

Relative Cathepsin B activity can be determined by comparing the Induced Sample relative fluorescence units (RFU) with the RFU values of the Uninduced Control or the Negative Control. If desired, the units of cathepsin B can be determined by generating a standard curve using free AFC (not included) under the assay conditions.

Frequently Asked Questions:

Can the kit work on bacteria or yeast cells?

The kit has been standardized for mammalian cells only.

Can an alternate buffer be used for sample preparation (cell lysis, sample dilutions, etc.)?

The Cathepsin B Activity Assay Kit assay buffers are optimized for the reactions the kit is designed for. The buffers not only contain detergents for efficient lysis of cells/tissue, but also contain proprietary components required for the further reactions. Therefore, it is highly recommended to use the buffers provided in the kit for the best results.



Notice

We provide information and advice to our customers on application technologies and regulatory matters to the best of our knowledge and ability, but without obligation or liability. Existing laws and regulations are to be observed in all cases by our customers. This also applies in respect to any rights of third parties. Our information and advice do not relieve our customers of their own responsibility for checking the suitability of our products for the envisaged purpose.

The information in this document is subject to change without notice and should not be construed as a commitment by the manufacturing or selling entity, or an affiliate. We assume no responsibility for any errors that may appear in this document.

Contact Information

For the location of the office nearest you, go to SigmaAldrich.com/offices.

Technical Service

Visit the tech service page on our web site at SigmaAldrich.com/techservice.

Standard Warranty

The applicable warranty for the products listed in this publication may be found at SigmaAldrich.com/terms.