



ab112154

Proteasome 20S Activity Assay Kit (Fluorometric)

Instructions for Use

For detecting proteasome activity in cultured cells using our proprietary green fluorescence probe.

This product is for research use only and is not intended for diagnostic use.



ab112154 Proteasome 20S Activity Assay Kit (Fluorometric)

Table of Contents

| | |
|-------------------------|----|
| 1. Introduction | 3 |
| 2. Protocol Summary | 5 |
| 3. Kit Contents | 6 |
| 4. Storage and Handling | 6 |
| 5. Assay Protocol | 7 |
| 6. Data Analysis | 10 |
| 7. Troubleshooting | 11 |

1. Introduction

The main function of the proteasome is to degrade unneeded or damaged proteins by proteolysis, a chemical reaction that breaks peptide bonds. The proteasomal degradation pathway is essential for many cellular processes, including the cell cycle, the regulation of gene expression, and the responses to oxidative stress. The most common form of the proteasome in this pathway is the proteasome 26S, an ATP-dependent proteolytic complex, which contains one 20S (700-kDa) core particle structure and two 19S (700-kDa) regulatory caps. The 20S core contains three major proteolytic activities including chymotrypsin-like, trypsin-like and caspase-like activities. It is responsible for the breakdown of the key proteins involved with apoptosis, DNA repair, endocytosis, and cell cycle control.

ab112154 Proteasome 20S Activity Assay Kit is a homogeneous fluorescent assay that measures the chymotrypsin-like protease activity associated with the proteasome complex in cultured cells. ab112154 uses LLVY-R110 as a fluorogenic indicator for proteasome activities. Cleavage of LLVY-R110 by the proteasome generates strongly green fluorescent R110 that is monitored fluorometrically at 520-530 nm with excitation at 480-500 nm. The kit provides all the essential components with an optimized assay protocol. The assay is robust, and can be readily adapted for high-

ab112154 Proteasome 20S Activity Assay Kit (Fluorometric)

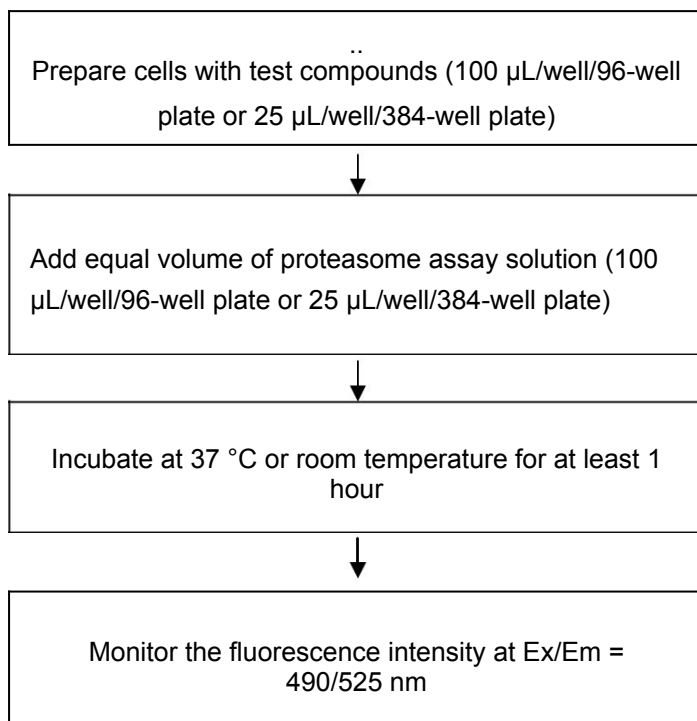
throughput assays to evaluate the proteasome activities or screen inhibitors in cultured cells or in solution. The assay can be performed in a convenient 96-well and 384-well fluorescence microtiter-plate format.

Kit Key Features

- Continuous:** Easily adapted to automation without a separation step.
- Convenient:** Includes all essential assay components.
- Increased Sensitivity:** Increased signal to background ratio.
- Versatile Applications:** Compatible with many cell lines and targets.

2. Protocol Summary

Summary for One 96-well Plate



Note: Thaw all the kit components to room temperature before starting the experiment.

3. Kit Contents

| Components | Amount |
|---|---------------|
| Component A: Proteasome LLVY-R110 Substrate | 1 vial |
| Component B: Assay Buffer | 1 x 10 mL |
| Component C: DMSO | 100 μ L |

4. Storage and Handling

Keep at -20 °C. Avoid exposure to light.

5. Assay Protocol

Note: *This protocol is for one 96 - well plate.*

A. Preparation of Cells

1. For adherent cells: Plate cells overnight in growth medium at 80,000 cells/well/90 μ L for a 96-well plate or 20,000cells/well/20 μ L for a 384-well plate.
2. For non-adherent cells: Centrifuge the cells from the culture medium and then suspend the cell pellet in culture medium at 300,000 cells/well/90 μ L for a 96-well poly-D lysine plate or 80,000 cells/well/20 μ L for a 384-well poly-D lysine plate. Centrifuge the plate at 800 rpm for 2 minutes with brake off prior to the experiments.

Note: Each cell line should be evaluated on an individual basis to determine the optimal cell density.

B. Preparation of Proteasome Assay Loading Solution

1. Thaw all the kit components at room temperature before use.
2. Make 400X Proteasome LLVY-R110 Substrate stock solution: Add 25 μ L of DMSO (Component C) to the vial

ab112154 Proteasome 20S Activity Assay Kit (Fluorometric)

of Proteasome LLVY-R110 Substrate (Component A), and mix well.

3. Make proteasome assay loading solution: Add 25 μL of 400X Proteasome LLVY-R110 Substrate stock solution (from Step 2) into 10 mL of Assay Buffer (Component B), and mix well.

Note: 25 μL of 400X Proteasome LLVY-R110 Substrate stock solution (from Step 2) and 10 mL of Assay Buffer (Component B) are enough for 1 plate. Aliquot and store the unused 400X Proteasome LLVY-R110 Substrate stock solution and Assay Buffer at $-20\text{ }^{\circ}\text{C}$. Avoid repeated freeze-thaw cycles.

C. Run Proteasome Assay:

1. Treat cells with 10 μL of 10X test compound (for a 96-well plate) or 5 μL of 5X test compound (for a 384-well plate) in PBS or desired buffer. For blank wells (medium without the cells), add the corresponding amount of compound buffer.
2. Incubate the cell plates in a 5% CO_2 , 37 $^{\circ}\text{C}$ incubator for a desired period of time.

ab112154 Proteasome 20S Activity Assay Kit (Fluorometric)

Note: Pure proteasome or cell lysates can be used directly for screening the proteasome inhibitors.

3. Add 100 μL /well (96-well plate) or 25 μL /well (384-well plate) of proteasome assay loading solution (from Step B.3).

4. Incubate the plate at 37 °C or room temperature for at least 1 hour (2 hours to overnight), protected from light.

Note: Each cell line should be evaluated on an individual basis to determine the optimal incubation time.

5. Monitor the fluorescence intensity (top read) at Ex/Em = 490/525 nm.

6. Data Analysis

The fluorescence in blank wells with the growth medium is subtracted from the values for those wells with the cells. The background fluorescence of the blank wells may vary depending on the sources of the growth media or the microtiter plates

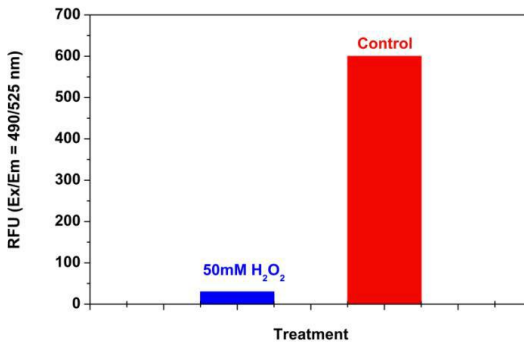


Figure 1. Detection of Proteasome Activity in Jurkat cells. Jurkat cells were seeded on the same day at 500,000 cells/90 μ L/well in a 96-well black wall/clear bottom plate. The cells were treated with or without 50 mM H₂O₂ for 30 minutes. The proteasome assay loading solution (100 μ L/well) was added and incubated in a 5% CO₂, 37 °C incubator for 3 hours. The fluorescence intensity was measured at Ex/Em = 490/525 using a fluorescent microplate reader.

7. Troubleshooting

| Problem | Reason | Solution |
|--------------------|--|---|
| Assay not working | Assay buffer at wrong temperature | Assay buffer must not be chilled - needs to be at RT |
| | Protocol step missed | Re-read and follow the protocol exactly |
| | Plate read at incorrect wavelength | Ensure you are using appropriate reader and filter settings (refer to datasheet) |
| | Unsuitable microtiter plate for assay | Fluorescence: Black plates (clear bottoms); Luminescence: White plates; Colorimetry: Clear plates. If critical, datasheet will indicate whether to use flat- or U-shaped wells |
| Unexpected results | Measured at wrong wavelength | Use appropriate reader and filter settings described in datasheet |
| | Samples contain impeding substances | Troubleshoot and also consider deproteinizing samples |
| | Unsuitable sample type | Use recommended samples types as listed on the datasheet |
| | Sample readings are outside linear range | Concentrate/ dilute samples to be in linear range |

ab112154 Proteasome 20S Activity Assay Kit (Fluorometric)

| | | |
|---|---|---|
| Samples with inconsistent readings | Unsuitable sample type | Refer to datasheet for details about incompatible samples |
| | Samples prepared in the wrong buffer | Use the assay buffer provided (or refer to datasheet for instructions) |
| | Samples not deproteinized (if indicated on datasheet) | Use the 10kDa spin column (ab93349) or Deproteinizing sample preparation kit (ab93299) |
| | Cell/ tissue samples not sufficiently homogenized | Increase sonication time/ number of strokes with the Dounce homogenizer |
| | Too many freeze-thaw cycles | Aliquot samples to reduce the number of freeze-thaw cycles |
| | Samples contain impeding substances | Troubleshoot and also consider deproteinizing samples |
| | Samples are too old or incorrectly stored | Use freshly made samples and store at recommended temperature until use |
| Lower/ Higher readings in samples and standards | Not fully thawed kit components | Wait for components to thaw completely and gently mix prior use |
| | Out-of-date kit or incorrectly stored reagents | Always check expiry date and store kit components as recommended on the datasheet |
| | Reagents sitting for extended periods on ice | Try to prepare a fresh reaction mix prior to each use |
| | Incorrect incubation time/ temperature | Refer to datasheet for recommended incubation time and/ or temperature |
| | Incorrect amounts used | Check pipette is calibrated correctly (always use smallest volume pipette that can pipette entire volume) |

ab112154 Proteasome 20S Activity Assay Kit (Fluorometric)

| Problem | Reason | Solution |
|------------------------------|--|--|
| Standard curve is not linear | Not fully thawed kit components | Wait for components to thaw completely and gently mix prior use |
| | Pipetting errors when setting up the standard curve | Try not to pipette too small volumes |
| | Incorrect pipetting when preparing the reaction mix | Always prepare a master mix |
| | Air bubbles in wells | Air bubbles will interfere with readings; try to avoid producing air bubbles and always remove bubbles prior to reading plates |
| | Concentration of standard stock incorrect | Recheck datasheet for recommended concentrations of standard stocks |
| | Errors in standard curve calculations | Refer to datasheet and re-check the calculations |
| | Use of other reagents than those provided with the kit | Use fresh components from the same kit |

For further technical questions please do not hesitate to contact us by email (technical@abcam.com) or phone (select “contact us” on www.abcam.com for the phone number for your region).

ab112154 Proteasome 20S Activity Assay Kit (Fluorometric)

Abcam in the USA

Abcam Inc
1 Kendall Square, Ste B2304
Cambridge,
MA 02139-1517
USA

Toll free: 888-77-ABCAM (22226)
Fax: 866-739-9884

Abcam in Europe

Abcam plc
330 Cambridge Science Park
Cambridge
CB4 0FL
UK

Tel: +44 (0)1223 696000
Fax: +44 (0)1223 771600

Abcam in Japan

Abcam KK
1-16-8 Nihonbashi
Kakigaracho,
Chuo-ku, Tokyo
103-0014
Japan

Tel: +81-(0)3-6231-094
Fax: +81-(0)3-6231-0941

Abcam in Hong Kong

Abcam (Hong Kong) Ltd
Unit 225A & 225B, 2/F
Core Building 2
1 Science Park West Avenue
Hong Kong Science Park
Hong Kong

Tel: Tel: 400 921 0189 / +86 21
2070 0500
Fax: (852) 3016-1888