

# Proteasome 20S Activity Assay Kit (Fluorometric) ab112154

[2 References](#) [画像数 1](#)

### 製品の概要

製品名	Proteasome 20S Activity Assay Kit (Fluorometric)
検出方法	Fluorescent
サンプルの種類	Adherent cells, Suspension cells
アッセイタイプ	Quantitative
種交差性	交差種: Mammals, Other species
製品の概要	<p>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-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.</p>

Visit our [FAQs page](#) for tips and troubleshooting.

### 特記事項

Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances. It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.

### 試験プラットフォーム

Microplate reader

### 製品の特性

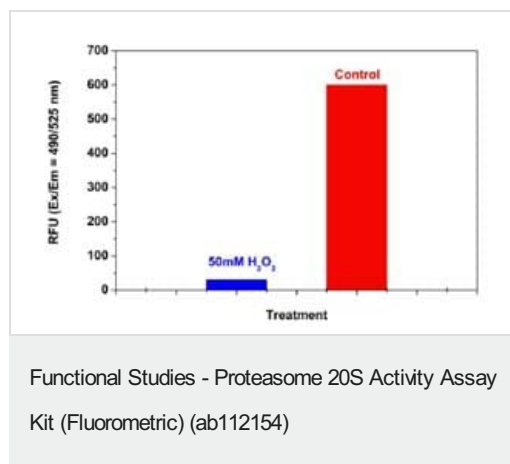
#### 保存方法

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

内容	ラベル	1 x 96 tests
Assay Buffer		1 x 10ml
DMSO		1 x 100µl

内容	ラベル	1 x 96 tests
Proteasome LLVY-R110 Substrate	Component A	1 vial

## 画像



Detection of Proteasome Activity in Jurkat cells. Jurkat cells were seeded on the same day at 500,000 cells/90  $\mu$ L/well in a 96-well black wall/clear bottom plate. The cells were treated with or without 50 mM H<sub>2</sub>O<sub>2</sub> for 30 minutes. The proteasome assay loading solution (100  $\mu$ L/well) was added and incubated in a 5% CO<sub>2</sub>, 37 °C incubator for 3 hours. The fluorescence intensity was measured at Ex/Em = 490/525.

**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.co.jp/abpromise> or contact our technical team.

## Terms and conditions

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