

# Proteasome Activity Assay Kit ab107921

★★★★★ [2 Abreviews](#) [69 References](#) [4 Images](#)

### Overview

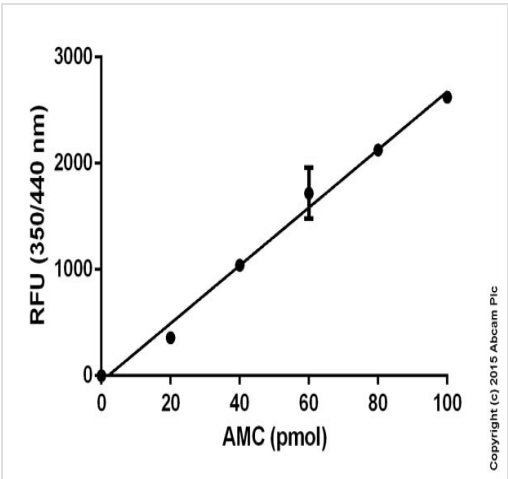
<b>Product name</b>	Proteasome Activity Assay Kit
<b>Detection method</b>	Fluorescent
<b>Sample type</b>	Cell Lysate
<b>Assay type</b>	Enzyme activity
<b>Assay time</b>	1h 00m
<b>Species reactivity</b>	<b>Reacts with:</b> Mammals, Other species
<b>Product overview</b>	<p>Proteasome Activity Assay Kit (ab107921) takes advantage of the chymotrypsin-like activity, using an AMC-tagged peptide substrate (Proteasome Substrate (Succ-LLVY-AMC in DMSO), which releases free, highly fluorescent AMC (Ex/Em 350/440 nm) in the presence of proteolytic activity.</p> <p>The kit also includes a positive control (Jurkat Cell lysate with significant proteasome activity) and a specific proteasome inhibitor MG-132 which suppresses all proteolytic activity due to proteasomes. This permits differentiation of proteasome activity from other protease activity which may be present in samples.</p> <p>Proteasome activity assay protocol summary:</p> <ul style="list-style-type: none"> <li>- add samples, standards and positive control to wells</li> <li>- add proteasome inhibitor to half of sample and positive control wells</li> <li>- add proteasome substrate to sample and control wells</li> <li>- incubate for 20 min</li> <li>- analyze with microplate reader, incubate for further 30 min, analyze again</li> </ul>
<b>Notes</b>	<p>This product is manufactured by BioVision, an Abcam company and was previously called K245 Proteasome Activity Fluorometric Assay Kit. K245-100 is the same size as the 100 test size of ab107921.</p> <p>The 20S proteasome assembly is the functional protease structure with chymotrypsin-like, trypsin-like and caspase-like protease activities</p> <p>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.</p> <p>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.</p>
<b>Platform</b>	Microplate reader

Properties

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

Components	100 tests
AMC Standard	1 x 100µl
Proteasome Assay Buffer	1 x 25ml
Proteasome Inhibitor	1 x 100µl
Proteasome Positive Control	1 vial
Proteasome Substrate	1 x 100µl

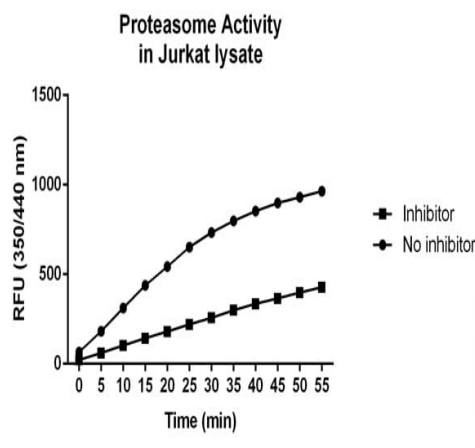
Images



Standard curve: mean of duplicates (+/- SD) with background reads subtracted

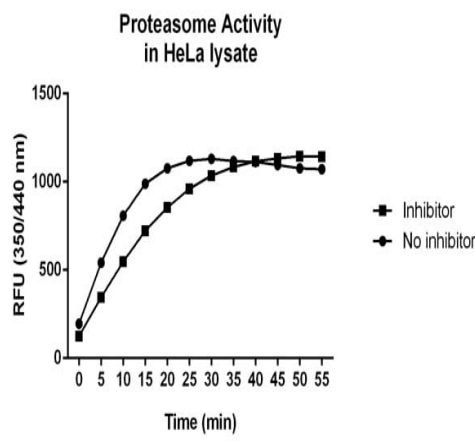
Functional Studies - Proteasome Activity Assay Kit  
(ab107921)

Proteasome Activity measured in Jurkat lysate



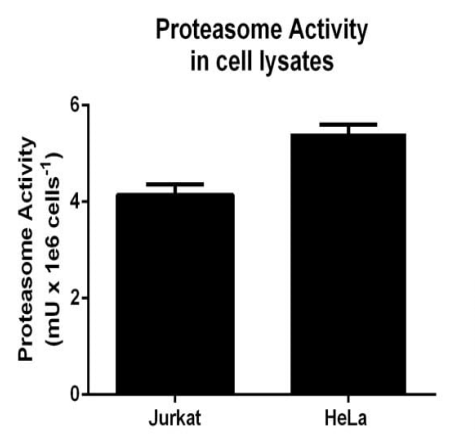
Functional Studies - Proteasome Activity Assay Kit  
(ab107921)

Proteasome Activity measured in HeLa lysate



Functional Studies - Proteasome Activity Assay Kit  
(ab107921)

Proteasome Activity measured in cell lysates showing activity (mU) per 1 mln cells.



Functional Studies - Proteasome Activity Assay Kit  
(ab107921)

Samples with the concentration of 3.3e6 cells/mL were used. 50 µl of each sample was used.

**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