

Proteasome 20S Activity Assay Kit (Fluorometric) ab112154

2 References 1 图像

概述

产品名称	Proteasome 20S Activity Assay试剂盒(Fluorometric)
检测方法	Fluorescent
样品类型	Adherent cells, Suspension cells
检测类型	Quantitative
种属反应性	与反应: Mammals, Other species
产品概述	<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.

说明	<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>
----	--

平台	Microplate reader
----	-------------------

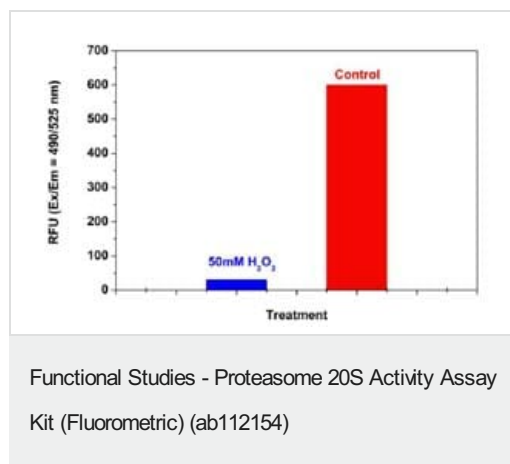
性能

存放说明	Store at -20°C. Please refer to protocols.
------	--

组件	标识符	1 x 96 tests
Assay Buffer		1 x 10ml
DMSO		1 x 100µl

组件	标识符	1 x 96 tests
Proteasome LLVY-R110 Substrate	Component A	1 vial

图片



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 H_2O_2 for 30 minutes. The proteasome assay loading solution (100 μ L/well) was added and incubated in a 5% CO_2 , 37 $^{\circ}C$ 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.cn/abpromise> or contact our technical team.

Terms and conditions

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