

Product datasheet

HIV-1 Protease Activity Assay Kit (Fluorometric)  
ab211105

4 Images

Overview

<b>Product name</b>	HIV-1 Protease Activity Assay Kit (Fluorometric)
<b>Detection method</b>	Fluorescent
<b>Sample type</b>	Purified protein, Tissue Lysate
<b>Assay type</b>	Quantitative
<b>Product overview</b>	HIV-1 Protease Activity Assay Kit (Fluorometric) (ab211105) provides a convenient and simple method for detecting the activity of recombinant and/or purified HIV-1 protease. The assay kit is based on the ability of active HIV-1 protease to cleave a synthetic peptide substrate to release the free fluorophore which can be easily quantified at Ex/Em = 330/450 nm using a fluorescence microplate reader.

**Notes** Human Immunodeficiency Virus (HIV) is the cause of the Acquired Immunodeficiency Syndrome (AIDS). HIV-1 protease is a retroviral aspartyl protease (retropepsin) that is essential for the life-cycle of the virus as it cleaves newly synthesized polyproteins at the appropriate places to create the mature protein components of an infectious HIV virion. Without effective HIV1 protease, HIV virions remain non-infectious.

**Platform** Microplate reader

Properties

**Storage instructions** Store at -80°C. Please refer to protocols.

Components	100 tests
Fl. Standard (10 mM in DMSO)	1 x 20µl
HIV-1 Protease (Positive Control)	1 x 7µl
HIV-1 Protease Assay Buffer	1 x 25ml
HIV-1 Protease Dilution Buffer	1 x 1ml

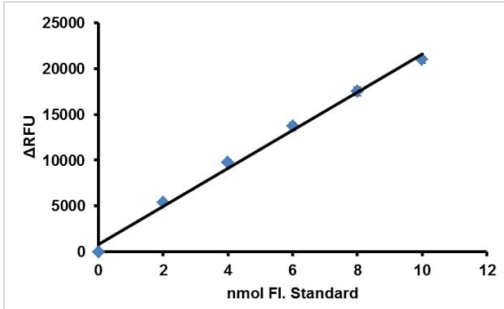
**Components**

**100 tests**

HIV-1 Protease Substrate

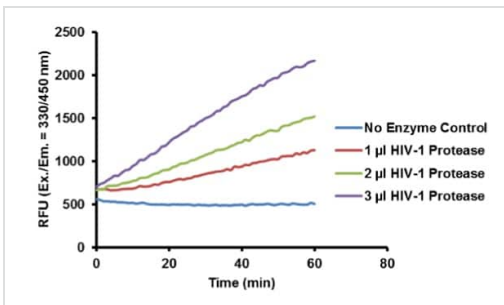
1 x 200µl

**Images**



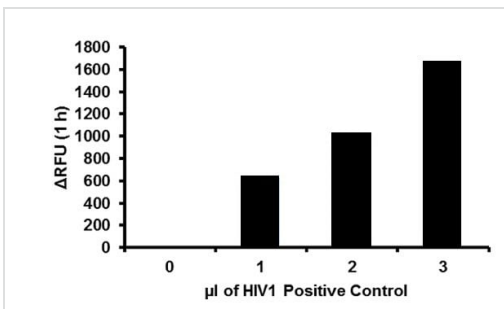
Typical FI. standard calibration curve.

HIV-1 Protease Activity Assay Kit (Fluorometric)  
(ab211105)



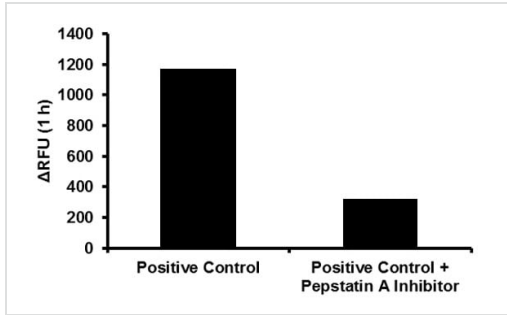
Kinetic progress curves observed for increasing amounts of HIV-1 Protease (positive control, included in the kit) through time.

HIV-1 Protease Activity Assay Kit (Fluorometric)  
(ab211105)



Measurement of HIV-1 Protease activity using increasing HIV-1 Protease positive control included in the kit, calculated from data from kinetic curves.

HIV-1 Protease Activity Assay Kit (Fluorometric)  
(ab211105)



Measurement of HIV-1 Protease activity using HIV-1 Protease included in kit as positive control and activity of HIV-1 Protease in presence of Pepstatin A (protease inhibitor).

HIV-1 Protease Activity Assay Kit (Fluorometric)  
(ab211105)

**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