

Product datasheet

HIV-1 Protease Inhibitor Assay Kit (Fluorometric)
ab211106

1 Image

Overview

Product name	HIV-1 Protease Inhibitor Assay Kit (Fluorometric)
Detection method	Fluorescent
Sample type	Inhibitor compounds
Assay type	Semi-quantitative
Product overview	HIV-1 Protease Inhibitor Screening Kit (Fluorometric) (ab211106) provides a rapid, simple test suitable for high-throughput screening of HIV-1 Protease inhibitors. The assay is based on the ability of active HIV-1 protease to cleave a synthetic peptide substrate to release a free fluorophore, which can be easily quantified using a fluorescence microplate reader. In the presence of a HIV-1 Protease inhibitor, the cleavage of the substrate is reduced/abolished, resulting in decrease or total loss of the fluorescence.

This simple and high-throughput adaptable assay kit can be used to screen/study/characterize potential inhibitors of HIV-1 Protease.

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 HIV-1 protease, HIV virions remain non-infectious.
--------------	---

Platform	Microplate reader
-----------------	-------------------

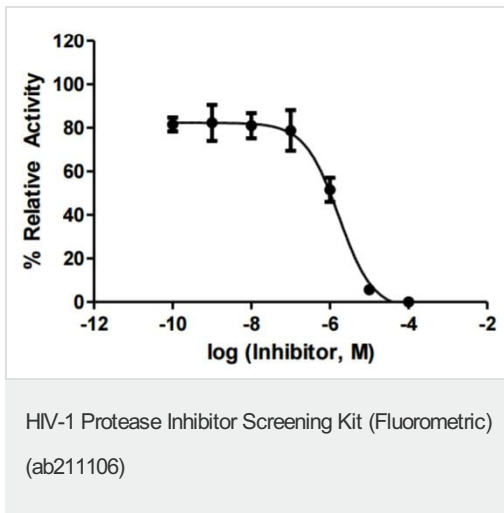
Properties

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

Components	100 tests
HIV-1 Protease	2 vials
HIV-1 Protease Assay Buffer	1 x 25ml
HIV-1 Protease Dilution Buffer	1 x 1ml

Components	100 tests
HIV-1 Protease Inhibitor (1 mM Pepstatin A in DMSO)	1 x 20µl
HIV-1 Protease Substrate	1 x 200µl

Images



Typical inhibition curve of HIV-1 Protease activity by the protease inhibitor Pepstatin A ($IC_{50} = 1.6 \mu M$) Assay was performed following the kit protocol.

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