

Product datasheet

Plasmin Inhibitor Screening Assay Kit (Fluorometric) ab204729

[2 Images](#)

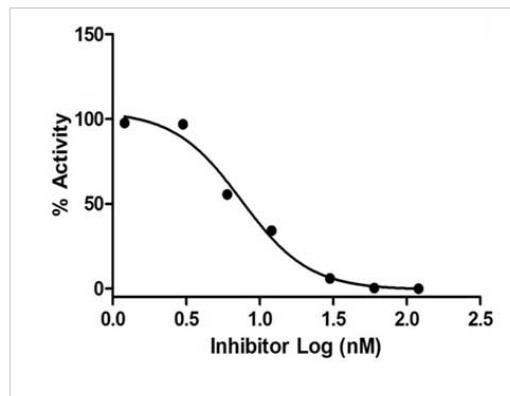
Overview

Product name	Plasmin Inhibitor Screening Assay Kit (Fluorometric)
Detection method	Fluorescent
Sample type	Inhibitor compounds
Assay type	Enzyme activity
Product overview	Plasmin Inhibitor Screening Assay Kit (Fluorometric) (ab204729) is based on the ability of Plasmin to cleave a synthetic AMC-based peptide substrate to release AMC, which can be detected by measuring its fluorescence at Ex/Em = 360/450 nm. In the presence of Plasmin specific inhibitors, the extent of cleavage reaction is reduced or completely abolished. The loss in the fluorescence intensity can be correlated to the amount of inhibitor present in the assay solution. The kit provides a simple and rapid method to screen potential inhibitors of Plasmin.
Notes	Plasmin (EC 3.4.21.7) is a serine protease occurring in plasma as plasminogen. Upon activation via cleavage by plasminogen activators; plasmin solubilizes fibrin clots and activates and/or degrades compounds of the coagulation and complement systems. Plasmin inhibitors are critical in the treatment of hyperfibrinolysis-associated blood loss and related complications.
Platform	Microplate reader

Properties

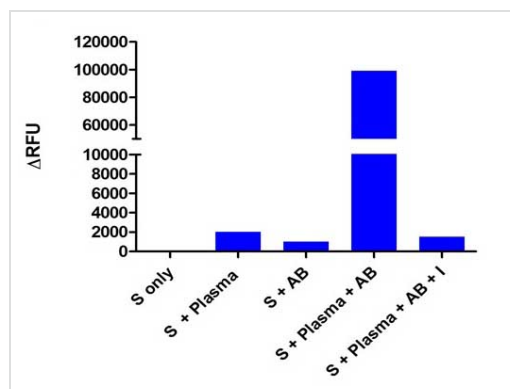
Storage instructions Store at -20°C. Please refer to protocols.

Components	100 tests
Plasmin Assay Buffer	1 x 15ml
Plasmin Dilution Buffer	1 x 1.5ml
Plasmin Enzyme	1 x 15µl
Plasmin Inhibitor (Aprotinin, 0.6 mM)	1 x 100µl
Plasmin Substrate	1 x 200µl



Inhibition of Plasmin activity by a Plasmin Inhibitor (Aprotinin).

Plasmin Inhibitor Screening Assay Kit (Fluorometric)
(ab204729)



Plasmin activity was measured in plasma samples in the presence and absence of a Plasmin inhibitor, Aprotinin. S = Substrate, I = Inhibitor, AB = Activation Buffer containing Urokinase.

Plasmin Inhibitor Screening Assay Kit (Fluorometric)
(ab204729)

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