

# Plasminogen activator Inhibitor Type 1 Human Chromogenic Activity Assay Kit ab108894

[1 Image](#)

### Overview

<b>Product name</b>	Plasminogen activator Inhibitor Type 1 Human Chromogenic Activity Assay Kit
<b>Detection method</b>	Colorimetric
<b>Sample type</b>	Cell culture supernatant, Serum, Plasma
<b>Assay type</b>	Quantitative
<b>Sensitivity</b>	0.57 AU/ml
<b>Range</b>	0.625 IU/ml - 20 IU/ml
<b>Product overview</b>	Plasminogen activator Inhibitor Type 1 Human Chromogenic Activity Kit (ab108894) has been developed to determine human type I plasminogen activator inhibitor (PAI-1) activity in cell culture supernatant, serum, and plasma samples.

A fixed amount of tissue-type plasminogen activator (tPA) is added in excess to undiluted sample, which allows PAI-1 and tPA to form an inactive complex. The assay measures plasminogen activation by residual tPA in coupled assays that contain tPA, plasminogen, and a plasmin-specific synthetic substrate. The amount of plasmin produced is quantitated using a highly specific plasmin substrate releasing a yellow para-nitroaniline (pNA) chromophore. The absorbance of the pNA at 405 nm is inversely proportional to the PAI-1 enzymatic activity.

**Upon arrival, immediately store components of the kit at recommended temperatures. Avoid repeated freeze-thaw cycles.**

<b>Notes</b>	Type I plasminogen activator inhibitor (PAI-1) is a 50 kDa serpin family member that inhibits tissue- and urokinase-type plasminogen activators (t-PA, u-PA). This protein appears to be an important regulator of plasminogen activation by t-PA and extracellular proteolysis by u-PA. The plasminogen activator proteolytic enzyme systems are important not only for fibrinolysis but also for extracellular matrix remodeling, and have been implicated in a number of normal and pathological processes including angiogenesis, ovulation and embryogenesis, thrombotic and hemorrhagic disorders, connective tissue diseases, neoplasm and sepsis. PAI-1 is a prognosticator in breast cancer, gastric cancer, various forms of lung cancer and cervical cancer.
--------------	---

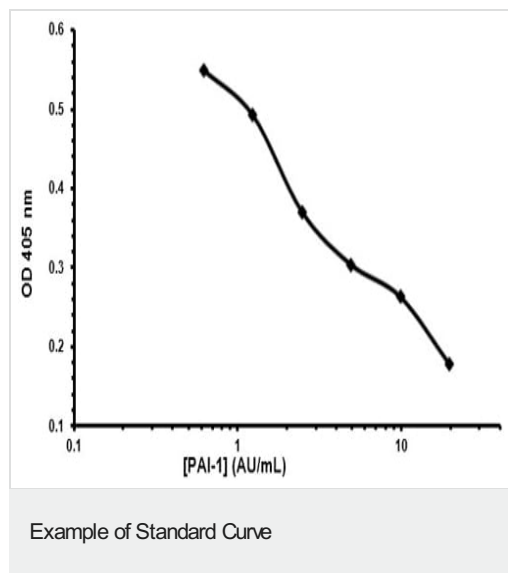
<b>Platform</b>	Microplate reader
-----------------	-------------------

### Properties

**Storage instructions**

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

Components	100 tests
Assay Diluent	1 x 30ml
Human Plasminogen (Lyophilized)	3 vials
Microplate	1 unit
Plasmin Substrate (Lyophilized)	2 vials
Sealing Tapes	3 units
tPA Standard 100IU (Lyophilized)	1 vial

**Images**

Representative Standard Curve using ab108894.

**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