

# Urokinase type Plasminogen Activator Human Chromogenic Activity Assay Kit - Direct Assay ab108916

1 [References](#) 2 [图像](#)

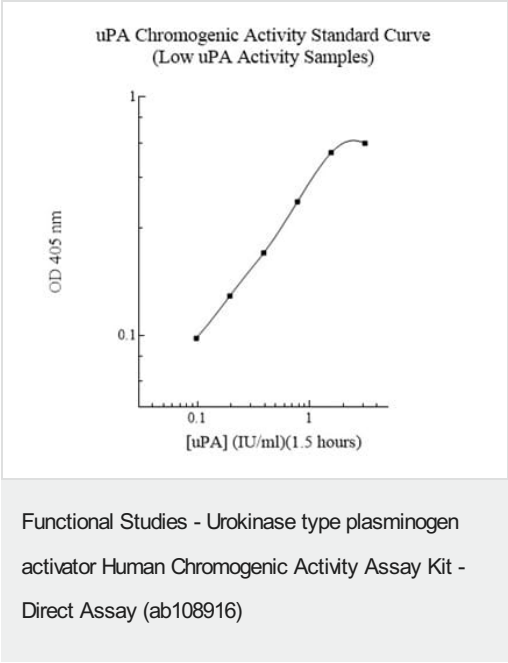
概述

产品名称	Urokinase type Plasminogen活化剂人Chromogenic Activity Assay试剂盒- Direct Assay
检测方法	Colorimetric
样品类型	Cell culture supernatant, Serum, Plasma
检测类型	Enzyme activity (quantitative)
灵敏度	< 0.062 IU/ml
种属反应性	与反应: Human
产品概述	<p>Urokinase type plasminogen activator (uPA) is a highly restricted serine protease that converts the zymogen plasminogen to active plasmin, a broad-spectrum serine proteinase capable of degrading most of the major protein components of the extracellular matrix. Binding of uPA to its receptor and subsequent uPA mediated pericellular proteolysis are involved in many process including cell migration and tissue remodeling in angiogenesis, atherogenesis, tumor cell metastasis, and ovulation. High level of uPA is a poor prognostic marker for aggressive breast cancer, aggressive prostate cancer, bladder cancer and gastric cancer.</p> <p>ab108916 uPA Human Chromogenic Activity Assay Kit is developed to determine Human uPA activity in plasma, serum and cell culture supernatants. The amidolytic activity of uPA is quantitated using a highly specific uPA substrate releasing a yellow para-nitroaniline (pNA) chromophore. The change in absorbance of the pNA at 405 nm is directly proportional to the uPA enzymatic activity.</p> <p><b>The entire kit may be stored at -20°C for long term storage before reconstitution - Avoid repeated freeze-thaw cycles.</b></p>
平台	Microplate reader
性能	
存放说明	Store at -20°C. Please refer to protocols.

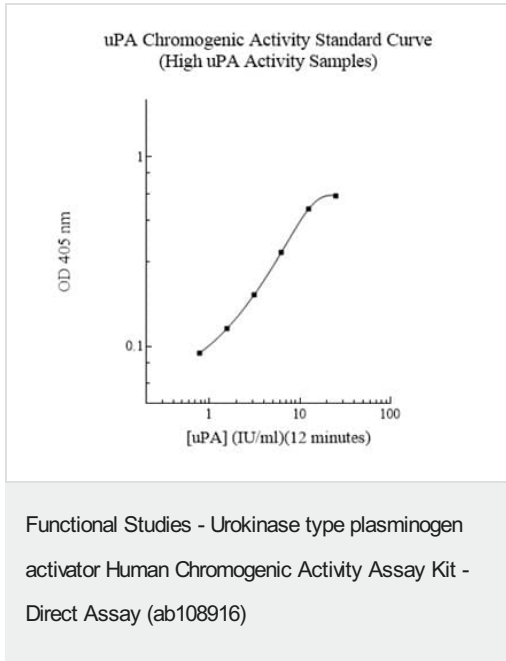
组件	96 tests
Assay Diluent	1 x 30ml
Human uPA Standard	1 vial
Microplate	1 unit
Sealing Tapes	3 units
uPA Substrate	3 units

功能	Specifically cleave the zymogen plasminogen to form the active enzyme plasmin.
组织特异性	Expressed in the prostate gland and prostate cancers.
序列相似性	<p>Belongs to the peptidase S1 family.</p> <p>Contains 1 EGF-like domain.</p> <p>Contains 1 kringle domain.</p> <p>Contains 1 peptidase S1 domain.</p>
翻译后修饰	Phosphorylation of Ser-158 and Ser-323 abolishes proadhesive ability but does not interfere with receptor binding.
细胞定位	Secreted.

图片



uPA Chromogenic Activity Standard Curve (For High uPA Activity Samples)



uPA Chromogenic Activity Standard Curve (For Low uPA Activity Samples)

**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