

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

[1 References](#) [画像数 2](#)

製品の概要

| | |
|------------|---|
| 製品名 | Urokinase type Plasminogen Activator Human Chromogenic Activity Assay Kit - Direct Assay |
| 検出方法 | Colorimetric |
| サンプルの種類 | Cell culture supernatant, Serum, Plasma |
| アッセイタイプ | Enzyme activity (quantitative) |
| 検出感度 | < 0.062 IU/ml |
| 種交差性 | 交差種: Human |
| 製品の概要 | <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>The entire kit may be stored at -20°C for long term storage before reconstitution - Avoid repeated freeze-thaw cycles.</p> |
| 試験プラットフォーム | Microplate reader |

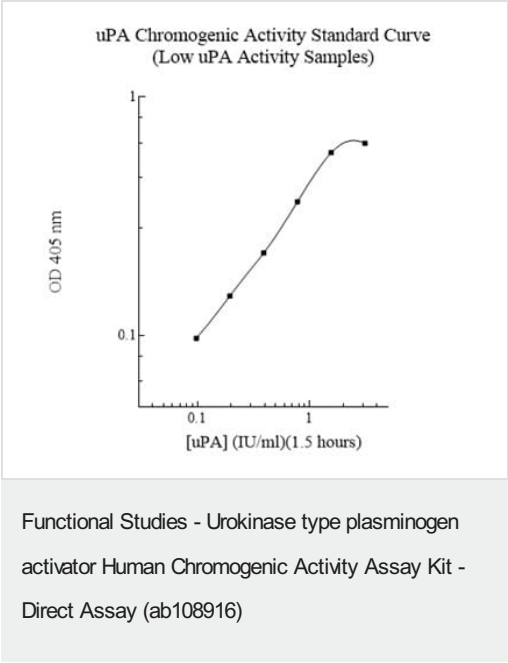
製品の特性

| | |
|------|--|
| 保存方法 | Store at -20°C. Please refer to protocols. |
|------|--|

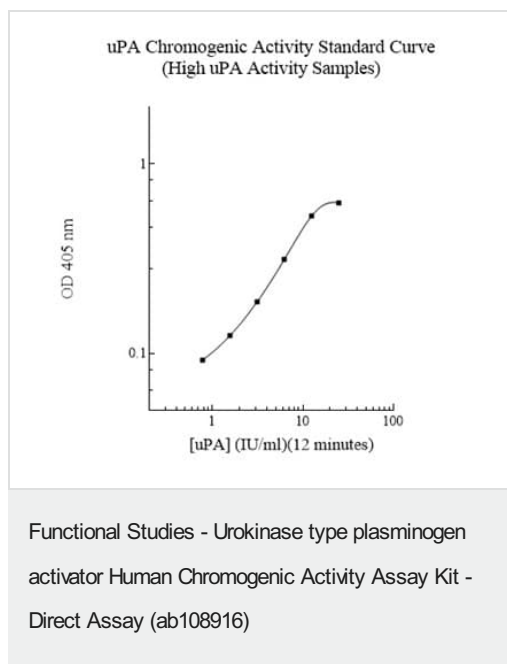
| 内容 | 96 tests |
|--------------------|----------|
| Assay Diluent | 1 x 30ml |
| Human uPA Standard | 1 vial |
| Microplate | 1 unit |
| Sealing Tapes | 3 units |
| uPA Substrate | 3 units |

| | |
|-------|--|
| 機能 | Specifically cleave the zymogen plasminogen to form the active enzyme plasmin. |
| 組織特異性 | Expressed in the prostate gland and prostate cancers. |
| 配列類似性 | <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> |
| 翻訳後修飾 | Phosphorylation of Ser-158 and Ser-323 abolishes proadhesive ability but does not interfere with receptor binding. |
| 細胞内局在 | Secreted. |

画像



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.co.jp/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors