

Urokinase Activity Assay Kit (Fluorometric) ab174098

[2 References](#) [3 Images](#)

Overview

Product name	Urokinase Activity Assay Kit (Fluorometric)
Detection method	Fluorescent
Sample type	Serum, Cell culture extracts, Nuclear Extracts, Purified protein
Assay type	Enzyme activity (quantitative)
Sensitivity	≥ 0.02 IU/ml
Species reactivity	Reacts with: Mammals, Other species
Product overview	Urokinase Activity Assay Kit (Fluorometric) (ab174098) uses an AMC-based peptide substrate containing the recognition sequence for Urokinase. Urokinase present in the sample catalyzes the cleavage of the substrate and releases AMC, which can be easily quantified by measuring its fluorescence at Ex/Em = 350/450 nm. This sensitive assay detects very low levels of Urokinase activity in biological samples as well as purified enzymes (as low as 0.02 IU/ml).

Visit our [FAQs](#) page for tips and troubleshooting.

Notes This product is manufactured by BioVision, an Abcam company and was previously called K728 Urokinase Activity Fluorometric Assay Kit. K728-100 is the same size as the 100 test size of ab174098.

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 remodelling 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.

Platform Microplate reader

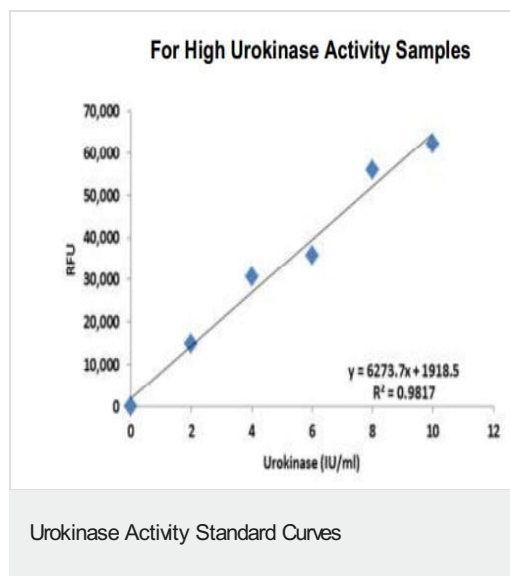
Properties

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

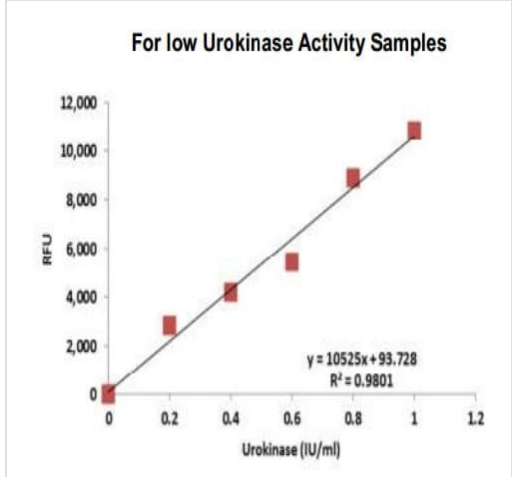
Components	
Cell Lysis Buffer	1 x 25ml
Human Urokinase (Standard/Positive Control)	1 x 1vial
Urokinase Assay Buffer	1 x 25ml
Urokinase Substrate	1 x 0.2ml

Function	Specifically cleave the zymogen plasminogen to form the active enzyme plasmin.
Tissue specificity	Expressed in the prostate gland and prostate cancers.
Sequence similarities	<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>
Post-translational modifications	Phosphorylation of Ser-158 and Ser-323 abolishes proadhesive ability but does not interfere with receptor binding.
Cellular localization	Secreted.

Images

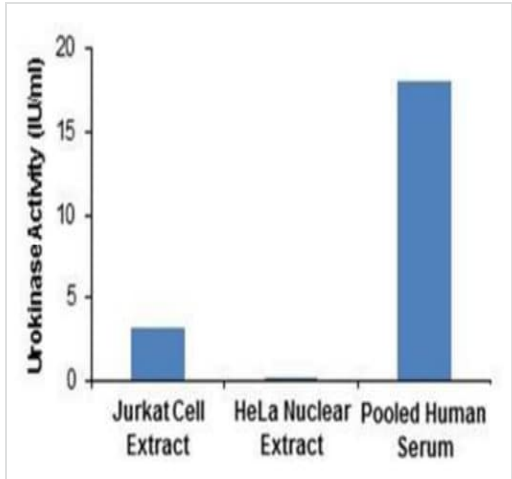


This is example data only.



This is example data only.

Urokinase Activity Standard Curves



Urokinase activities of Jurkat cell extract [10 µl], Hela nuclear extract [10 µl] and Pooled human serum (10 µl). Assays were performed following the kit protocol. This is example data only.

Urokinase activities in various 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.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