

Human uPA Receptor/U-PAR ELISA Kit (URKR) ab119612

2 References 1 Image

Overview

Product name	Human uPA Receptor/U-PAR ELISA Kit (URKR)
Detection method	Colorimetric
Sample type	Cell culture supernatant, Serum, Plasma, Other biological fluids, Tissue Extracts
Assay type	Sandwich (quantitative)
Sensitivity	< 4 pg/ml
Range	62.5 pg/ml - 4000 pg/ml
Assay duration	Multiple steps standard assay
Species reactivity	Reacts with: Human
Product overview	Abcam's Human uPA Receptor/U-PAR (URKR) <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate quantitative measurement of Human uPA Receptor/U-PAR (URKR) in cell culture supernatants, serum, plasma (heparin, EDTA) and urine.

A uPA Receptor/U-PAR (URKR) specific mouse monoclonal antibody has been precoated onto 96-well plates. Standards and test samples are added to the wells and incubated. A biotinylated detection polyclonal antibody from goat specific for uPA Receptor/U-PAR(URKR) is then added followed by washing with PBS or TBS buffer. Avidin-Biotin-Peroxidase Complex is added and unbound conjugates are washed away with PBS or TBS buffer. TMB is then used to visualize the HRP enzymatic reaction. TMB is catalyzed by HRP to produce a blue color product that changes into yellow after adding acidic stop solution. The density of yellow coloration is directly proportional to the Human uPA Receptor amount of sample captured in plate.

PlatformMicroplate

Properties

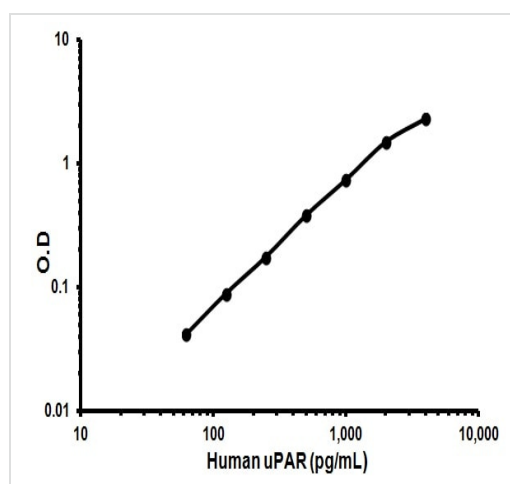
Storage instructionsStore at -20°C. Please refer to protocols.

Components	Identifier	1 x 96 tests
ABC Diluent Buffer	Blue Cap	1 x 12ml
Antibody Diluent Buffer	Green Cap	1 x 12ml

Components	Identifier	1 x 96 tests
Anti-Human uPA Receptor antibody Microplate (12 x 8 wells)		1 unit
Avidin-Biotin-Peroxidase Complex (ABC)		1 x 100µl
Biotinylated anti-Human uPA Receptor antibody		1 x 100µl
Lyophilized recombinant Human uPA Receptor standard		2 x 10ng
Plate Seal		1 x 4 units
Sample Diluent Buffer	Green Cap	1 x 30ml
TMB Color Developing Agent	Amber Bottle	1 x 10ml
TMB Stop Solution	Yellow Cap	1 x 10ml
Wash Buffer (25X)		1 x 20ml

Function	Acts as a receptor for urokinase plasminogen activator. Plays a role in localizing and promoting plasmin formation. Mediates the proteolysis-independent signal transduction activation effects of U-PA. It is subject to negative-feedback regulation by U-PA which cleaves it into an inactive form.
Tissue specificity	Expressed in neurons of the rolandic area of the brain (at protein level). Expressed in the brain.
Sequence similarities	Contains 3 UPAR/Ly6 domains.
Cellular localization	Secreted and Cell membrane.

Images



Typical Standard Curve

Representative standard curve using ab119612

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