

Human TAFI ELISA kit ab272774

1 Image

Overview

| | |
|--------------------|--|
| Product name | Human TAFI ELISA kit |
| Detection method | Colorimetric |
| Sample type | Serum, EDTA Plasma, Cit plasma |
| Assay type | Quantitative |
| Range | 1 ng/ml - 500 ng/ml |
| Assay duration | Multiple steps standard assay |
| Species reactivity | Reacts with: Human |
| Product overview | <p>Human TAFI ELISA kit (ab272774) is intended for the quantitative determination of total TAFI antigen in human plasma and serum.</p> <p>Human TAFI will bind to the affinity purified capture antibody coated on the microtiter plate. TAFI and TAFIa will react with the antibody on the plate. After appropriate washing steps, biotinylated primary antibody binds to the captured protein. Excess primary antibody is washed away and bound antibody is reacted with horseradish peroxidase-conjugated streptavidin. TMB substrate is used for color development at 450nm. A standard calibration curve is prepared along with the samples to be measured using dilutions of human TAFI. Color development is directly proportional to the concentration of total TAFI in the samples.</p> |
| Platform | Microplate |

Properties

Storage instructionsStore at +4°C. Please refer to protocols.

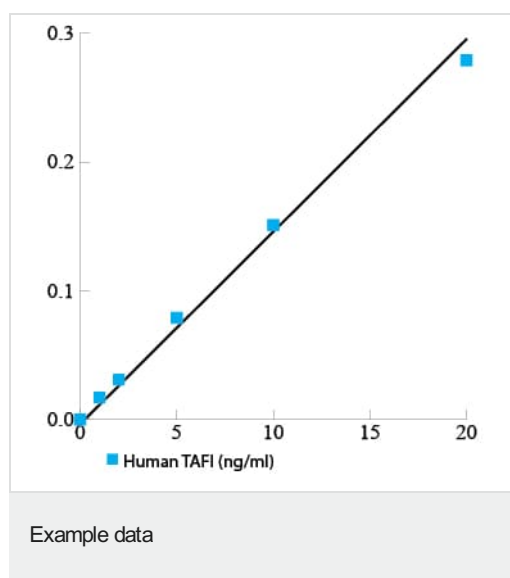
| Components | 1 x 96 tests |
|---|--------------|
| 10X Wash Buffer | 1 x 50ml |
| Anti-Human TAFI Primary Antibody Lyophilized Vial | 1 vial |
| HRP-conjugated Streptavidin | 1 vial |
| Human TAFI Standard Lyophilized Vial | 1 vial |
| | |

| Components | 1 x 96 tests |
|------------------|--------------|
| TAFI ELISA Plate | 1 unit |
| TMB Substrate | 1 x 10ml |

Relevance The thrombin-activatable fibrinolysis inhibitor (TAFI) is a hepatically secreted zymogen. It potently attenuates fibrinolysis by removing the fibrin C-terminal lysine and arginine residues that are important for the binding and activation of plasminogen. This molecule is also known as carboxypeptidase U (CPU). It is an unstable basic carboxypeptidase that circulates in human plasma in its proenzyme form.

Cellular localization Secreted

Images



A typical standard curve. Example only.

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