

Product datasheet

Human ANGPT4 ELISA Kit ab193693

[2 Images](#)

Overview

Product name Human ANGPT4 ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
--------	---	------	----	-----

Overall < 10%

Inter-assay

Sample	n	Mean	SD	CV%
--------	---	------	----	-----

Overall < 12%

Sample type Cell culture supernatant, Serum, Plasma

Assay type Sandwich (quantitative)

Sensitivity 0.25 ng/ml

Range 0.25 ng/ml - 200 ng/ml

Recovery

Sample specific recovery

Sample type	Average %	Range
Serum	101.4	88% - 120%
Plasma	83.13	67% - 91%
Cell culture media	113.7	95% - 130%

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview Abcam's ANGPT4 Human ELISA Kit (ab193693) is an in vitro enzyme-linked immunosorbent assay for the quantitative measurement of human ANGPT4 in serum plasma and cell culture supernatants.

This assay employs an antibody specific for human ANGPT4 coated on a 96-well plate. Standards and samples are pipetted into the wells and the immobilized antibody captures ANGPT4 present in the samples. The wells are washed and biotinylated anti-human ANGPT4 antibody is added. After washing away any unbound biotinylated antibody, an HRP-conjugated streptavidin is pipetted to the wells. After incubation, the wells are again washed, followed by the addition of a TMB substrate solution to the wells. Color will develop in proportion to the amount of ANGPT4 bound in each well. Addition of the Stop Solution will change the color from blue to yellow, and the intensity of the color is measured at 450 nm.

Platform Pre-coated microplate (12 x 8 well strips)

Properties

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

Components	1 x 96 tests
20X Wash Buffer Concentrate	1 x 25ml
5X Assay Diluent B	1 x 15ml
600X HRP-Streptavidin Concentrate	1 x 200µl
Assay Diluent A	1 x 30ml
Biotinylated Human Angiotensin-4 detection antibody (lyophilized)	2 vials
Human Angiotensin-4 Standards (lyophilized)	2 vials
Pre-coated Human Angiotensin-4 Microplate (12 strips x 8 wells)	1 unit
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

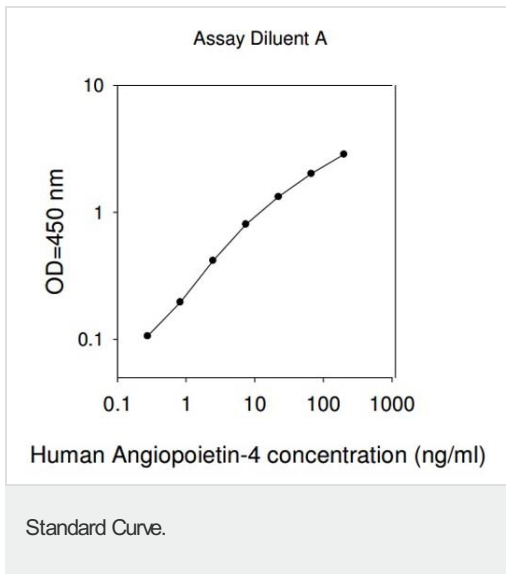
Function Binds to TEK/TIE2, modulating ANGPT1 signaling. Can induce tyrosine phosphorylation of TEK/TIE2. Promotes endothelial cell survival, migration and angiogenesis.

Tissue specificity Highly expressed in the lung with much lower levels found in other tissues.

Sequence similarities Contains 1 fibrinogen C-terminal domain.

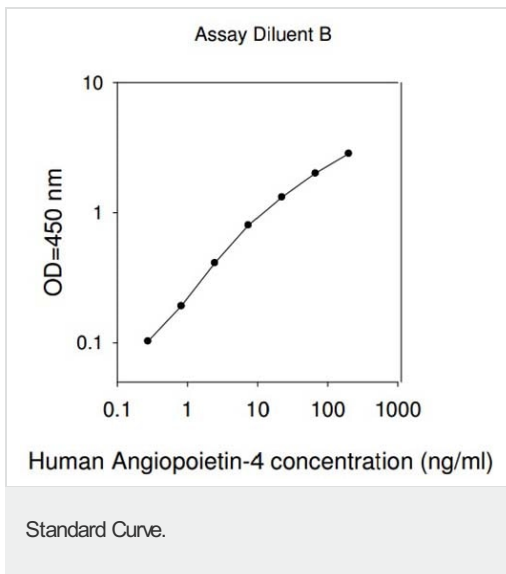
Cellular localization Secreted.

Images



Standard Curve in Assay Diluent A using ab193693 ANGPT4 Human ELISA Kit.

Data provided for demonstration purposes only. A new standard curve must be generated for each assay performed.



Standard Curve in Assay Diluent B using ab193693 ANGPT4 Human ELISA Kit.

Data provided for demonstration purposes only. A new standard curve must be generated for each assay performed.

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