

Human Angiopoietin 1 ELISA Kit (ANG1) ab99972

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

Overview

Product name	Human Angiopoietin 1 ELISA Kit (ANG1)
Detection method	Colorimetric
Sample type	Cell culture supernatant, Serum, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	< 30 pg/ml
Range	74 pg/ml - 18000 pg/ml
Recovery	< 95 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	95.25	83% - 105%
Serum	94.47	81% - 103%
Plasma	93.93	80% - 102%

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview Abcam's Angiopoietin 1 (ANG1) Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Human Angiopoietin 1 in serum, plasma, and cell culture supernatants.

This assay employs an antibody specific for Human Angiopoietin 1 coated on a 96-well plate. Standards and samples are pipetted into the wells and Angiopoietin 1 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human Angiopoietin 1 antibody is added. After washing away unbound biotinylated antibody, HRP-conjugated streptavidin is pipetted to the wells. The wells are again washed, a TMB substrate solution is added to the wells and color develops in proportion to the amount of Angiopoietin 1 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

Notes Optimisation may be required with urine samples.

Platform Microplate

Properties

Storage instructions

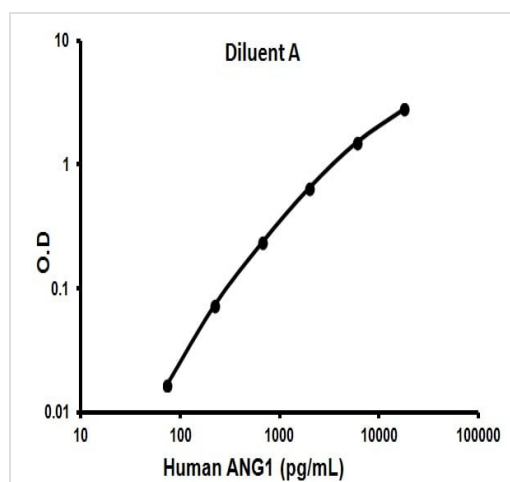
Store at -20°C. Please refer to protocols.

Components	1 x 96 tests
20X Wash Buffer	1 x 25ml
400X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent B	1 x 15ml
Angiopoietin 1 Microplate (12 x 8 well strips)	1 unit
Assay Diluent A	1 x 30ml
Biotinylated anti-Human Angiopoietin 1	2 vials
Recombinant Angiopoietin 1 Standard Human (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

Relevance

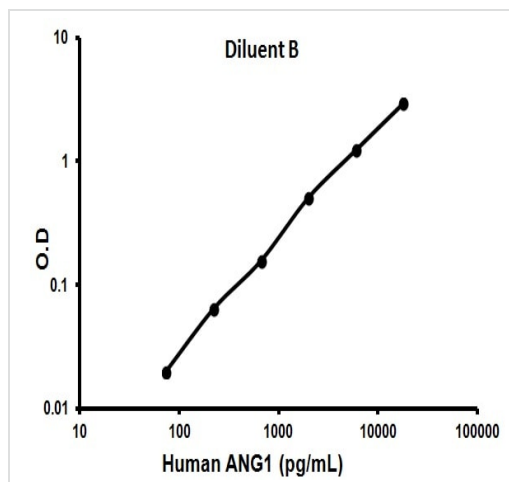
Angiopoietin 1 is one of the Angiopoietin (Ang) family of growth factors. It binds and activates the TIE2 receptor tyrosine kinase by inducing its tyrosine phosphorylation. It is implicated in endothelial developmental processes later and distinct from that of VEGF. The structure of angiopoietin 1 differs from that of known angiogenic factors or other ligands for receptor tyrosine kinases. Angiopoietin 1 appears to play a crucial role in mediating reciprocal interactions between the endothelium and surrounding matrix and mesenchyme. It also mediates blood vessel maturation/stability and it may play an important role in the heart early development. Angiopoietin 1 may have a potential therapeutic utility since it can be used for specifically targeting tumor vasculature or for promoting angiogenic processes in certain organs such as an ischemic heart.

Images



TYPICAL STANDARD CURVE - Assay Diluent A – Data provided for demonstration purposes only. A new standard curve must be generated for each assay performed.

TYPICAL STANDARD CURVE - Assay Diluent A



TYPICAL STANDARD CURVE - Assay Diluent B – Data provided for demonstration purposes only. A new standard curve must be generated for each assay performed.

TYPICAL STANDARD CURVE - Assay Diluent B

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