

Human Angiopoietin 2 ELISA Kit (ANG2) ab99971

[3 References](#) [1 Image](#)

Overview

Product name	Human Angiopoietin 2 ELISA Kit (ANG2)
Detection method	Colorimetric
Sample type	Cell culture supernatant, Serum, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	< 10 pg/ml
Range	4.12 pg/ml - 3000 pg/ml
Recovery	88 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	82.1	76% - 92%
Serum	85.3	82% - 109%
Plasma	96.9	95% - 116%

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview Abcam's Angiopoietin 2 (ANG2) Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Human Angiopoietin 2 in serum, plasma and cell culture supernates.

This assay employs an antibody specific for Human Angiopoietin 2 coated on a 96-well plate. Standards and samples are pipetted into the wells and Angiopoietin 2 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human Angiopoietin 2 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 2 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
200X HRP-Streptavidin Concentrate	1 x 200µl
20X Wash Buffer	1 x 25ml
5X Assay Diluent	1 x 15ml
Angiopoietin 2 Microplate (12 x 8 well strip)	1 unit
Biotinylated anti-Human Angiopoietin 2	2 vials
Recombinant Angiopoietin 2 Standard Human (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

Function

Can induce tyrosine phosphorylation of TIE2. Binds to TIE2 receptor and counteracts blood vessel maturation/stability mediated by angiopoietin-1. Its function may be context-dependent. In the absence of angiogenic inducers, such as VEGF, ANG2-mediated loosening of cell-matrix contacts may induce endothelial cell apoptosis with consequent vascular regression. In concert with VEGF, it may facilitate endothelial cell migration and proliferation, thus serving as a permissive angiogenic signal.

Sequence similarities

Contains 1 fibrinogen C-terminal domain.

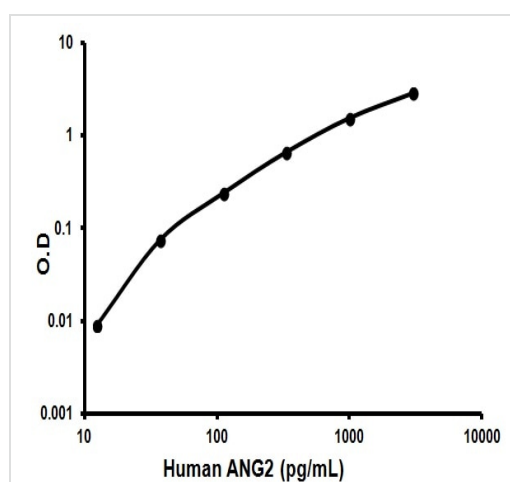
Domain

The Fibrinogen C-terminal domain mediates interaction with the TEK/TIE2 receptor.

Cellular localization

Secreted.

Images



Representative Standard Curve using ab99971.

Typical Standard Curve

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