

Human Angiopoietin 2 ELISA Kit (ANG2) ab99971

3 References 1 图像

概述

| | |
|------|---|
| 产品名称 | 人Angiopoietin 2 ELISA试剂盒(ANG2) |
| 检测方法 | Colorimetric |
| 样品类型 | Cell culture supernatant, Serum, Plasma |
| 检测类型 | Sandwich (quantitative) |
| 灵敏度 | < 10 pg/ml |
| 范围 | 4.12 pg/ml - 3000 pg/ml |
| 回收率 | 88 % |

特定样本回收率

| 样品类型 | 平均% | 范围 |
|--------------------------|------|------------|
| Cell culture supernatant | 82.1 | 76% - 92% |
| Serum | 85.3 | 82% - 109% |
| Plasma | 96.9 | 95% - 116% |

| | |
|-------|---|
| 实验步骤 | Multiple steps standard assay |
| 种属反应性 | 与反应: Human |
| 产品概述 | <p>Abcam's Angiopoietin 2 (ANG2) Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an <i>in vitro</i> enzyme-linked immunosorbent assay for the quantitative measurement of Human Angiopoietin 2 in serum, plasma and cell culture supernates.</p> <p>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.</p> |
| 说明 | Optimisation may be required with urine samples. |
| 平台 | Microplate |

性能

存放说明 Store at -20°C. Please refer to protocols.

| 组件 | 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 |

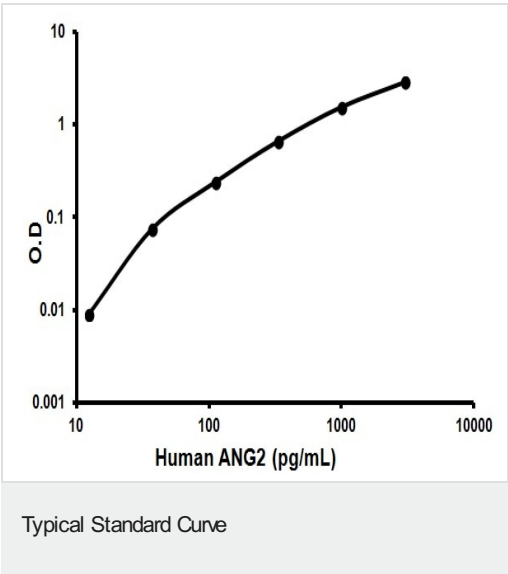
功能 Can induce tyrosine phosphorylation of TIE2. Binds to TIE2 receptor and counteracts blood vessel maturation/stability mediated by angiopoietin-1. Its function may be context-dependent. In the absence of angiogenic inducers, such as VEGF, ANG2-mediated loosening of cell-matrix contacts may induce endothelial 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.

序列相似性 Contains 1 fibrinogen C-terminal domain.

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

细胞定位 Secreted.

图片



Representative Standard Curve using ab99971.

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.cn/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors