

## Product datasheet

# Human Angiogenin ELISA Kit (ANG) ab99970

2 References 2 图像

### 概述

**产品名称** 人Angiogenin ELISA试剂盒(ANG)

**检测方法** Colorimetric

**精确度**

批次内				
样品	n	Mean	SD	CV%
Overall				< 10%

批次间				
样品	n	Mean	SD	CV%
Overall				< 12%

**样品类型** Cell culture supernatant, Serum, Plasma

**检测类型** Sandwich (quantitative)

**灵敏度** < 1.5 pg/ml

**范围** 1.64 pg/ml - 400 pg/ml

**复检** 94.7 %

特定样本复检结果		
样品类型	平均%	范围
Cell culture supernatant	95.37	85% - 104%
Serum	95.95	82% - 105%
Plasma	93.49	83% - 104%

**实验步骤** Multiple steps standard assay

**种属反应性** 与反应: Human

**产品概述** Abcam's Angiogenin (ANG) Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Human Angiogenin in serum, plasma and cell culture supernates.

This assay employs an antibody specific for Human Angiogenin coated on a 96-well plate. Standards and samples are pipetted into the wells and Angiogenin present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human Angiogenin 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 Angiogenin bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

**说明** Optimisation may be required with urine samples.

**经测试应用** **适用于:** Sandwich ELISA

**平台** Microplate

## 性能

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

组件	1 x 96 tests
20X Wash Buffer Concentrate	1 x 25ml
5X Assay Diluent B	1 x 15ml
700X HRP-Streptavidin Concentrate	1 x 200µl
Angiogenin Microplate (12 x 8 well strip)	1 unit
Assay Diluent A	1 x 30ml
Biotinylated anti-Human Angiogenin (lyophilized)	2 vials
Recombinant Human Angiogenin Standard	2 vials
Stop Solution	1 x 8ml
TMB Substrate Reagent	1 x 12ml

**相关性** Angiogenin is a member of the ribonuclease superfamily with approximately 35% amino acid sequence identity with pancreatic RNase. Like other members of the ribonuclease superfamily, angiogenin is a cytotoxic agent that can eliminate cellular protein synthesis. Angiogenin causes inhibition of protein synthesis by functioning as a cytotoxic tRNA-specific RNase. Angiogenesis is the formation of blood vessels or capillaries from existing blood vessels due to specific signals. Angiogenin induces the growth of new blood vessels. It stimulates capillary and umbilical vein endothelial cells to produce diacylglycerol and secrete prostacyclin by phospholipase activation. Angiogenin exhibits both angiogenic and non-angiogenic activities and is involved in endothelial cell migration, proliferation, and differentiation. It is produced by a variety of tumor and normal cell types. Cells that express angiogenin include vascular endothelial and smooth muscle cells, fibroblasts, normal colonic epithelium, normal peripheral blood lymphocytes, lung and colonic epithelial tumor cell lines, and primary gastrointestinal adenocarcinomas. Angiogenin is also present in normal human plasma.

**细胞定位** Nuclear and Secreted

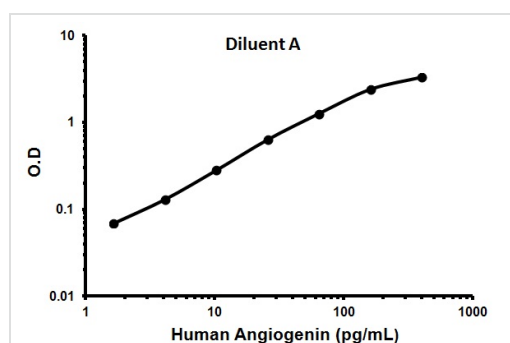
## 应用

Our [Abpromise guarantee](#) covers the use of **ab99970** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

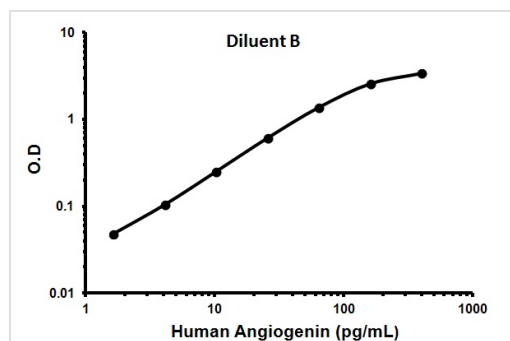
应用	Ab评论	说明
Sandwich ELISA		Use at an assay dependent concentration.

## 图片



Typical Standard Curve

Representative standard curve using ab99970.



Typical Standard Curve

Representative standard curve using ab99970.

**Please note:** All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

## 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 <http://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