

Human EG-VEGF ELISA Kit (PK1) ab119591

1 图像

概述

产品名称	人EG-VEGF ELISA试剂盒(PK1)
检测方法	Colorimetric
样品类型	Cell culture supernatant, Plasma
检测类型	Sandwich (quantitative)
灵敏度	< 7 pg/ml
范围	15.6 pg/ml - 1000 pg/ml
实验步骤	Multiple steps standard assay
种属反应性	与反应: Human
产品概述	Abcam's Human EG-VEGF (PK1) <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate quantitative measurement of Human EG-VEGF (PK1) in cell culture supernatants and plasma (heparin, EDTA).

A EG-VEGF (PK1) specific mouse monoclonal antibody has been precoated onto 96-well plates. Standards and test samples are added to the wells and incubated. A biotinylated detection polyclonal antibody from goat, specific for EG-VEGF (PK1) is then added followed by washing with PBS or TBS buffer. Avidin-Biotin-Peroxidase Complex is added and unbound conjugates are washed away with PBS or TBS buffer. TMB is then used to visualize the HRP enzymatic reaction. TMB is catalyzed by HRP to produce a blue color product that changes into yellow after adding acidic stop solution. The density of yellow coloration is directly proportional to the Human EG-VEGF (PK1) amount of sample captured in plate.

说明	This kit may not be sensitive enough to detect EG-VEGF in serum samples and this sample type is not covered by our AbPromise.
平台	Microplate

性能

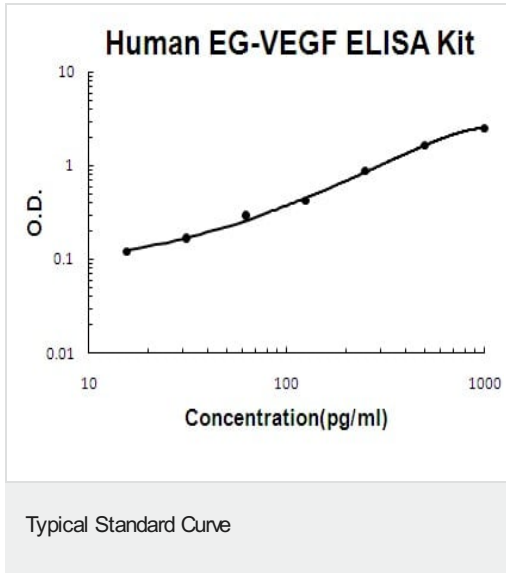
存放说明 Store at -20°C. Please refer to protocols.

组件	标识符	1 x 96 tests
ABC Diluent Buffer	Blue Cap	1 x 12ml

组件	标识符	1 x 96 tests
Antibody Diluent Buffer	Green Cap	1 x 12ml
Anti-Human EG-VEGF antibody Microplate (12 x 8 wells)		1 unit
Avidin-Biotin-Peroxidase Complex (ABC)		1 x 100µl
Biotinylated anti- Human EG-VEGF antibody		1 x 100µl
Lyophilized recombinant Human EG-VEGF standard		2 x 10ng
Plate Seal		1 x 4 units
Sample Diluent Buffer	Green Cap	1 x 30ml
TMB Color Developing Agent	Amber Bottle	1 x 10ml
TMB Stop Solution	Yellow Cap	1 x 10ml
Wash Buffer (25X)		1 x 20ml

功能	Potently contracts gastrointestinal (GI) smooth muscle. Induces proliferation, migration and fenestration (the formation of membrane discontinuities) in capillary endothelial cells derived from endocrine glands. Has little or no effect on a variety of other endothelial and non-endothelial cell types. Induces proliferation and differentiation, but not migration, of enteric neural crest cells. Directly influences neuroblastoma progression by promoting the proliferation and migration of neuroblastoma cells. Positively regulates PTGS2 expression and prostaglandin synthesis. May play a role in placentation. May play a role in normal and pathological testis angiogenesis.
组织特异性	Localizes to glandular epithelium, stroma and vascular epithelial cells of first trimester decidua (at protein level). Up-regulated in first trimester decidua when compared with non-pregnant endometrium. Expressed in the steroidogenic glands, ovary, testis, adrenal and placenta.
序列相似性	Belongs to the AVIT (prokineticin) family.
发展阶段	In adult testis, is strongly expressed only in Leydig cells. In testicular tumors, expressed specifically in Leydig cell tumors (at protein level). Expressed from 14 weeks until birth in fetal testis.
细胞定位	Secreted.

图片



Representative Standard Curve using ab119591.

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