

Human VEGF R1 ELISA Kit (FLT1) ab119613

3 References 1 图像

概述

产品名称 人VEGF R1 ELISA试剂盒(FLT1)

检测方法 Colorimetric

精确度 批次内

样品	n	Mean	SD	CV%
1	16	237pg/ml	9.71	= 4.1%
2	16	1369pg/ml	106.78	= 7.8%
3	16	4420pg/ml	203.32	= 4.6%

批次间

样品	n	Mean	SD	CV%
1	24	240pg/ml	11.28	= 4.7%
2	24	1352pg/ml	131.14	= 9.7%
3	24	4316pg/ml	202.85	= 4.7%

样品类型 Cell culture supernatant, Serum, Plasma

检测类型 Sandwich (quantitative)

灵敏度 < 30 pg/ml

范围 156 pg/ml - 10000 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 soluble VEGF R1 (FLT1) *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate quantitative measurement of Human soluble VEGF R1 (FLT1) in cell culture supernatants, serum and plasma (EDTA).

A VEGF R1 (FLT1) 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 VEGF R1 (FLT1) 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 VEGF R1 (FLT1) amount of sample captured in plate.

Get higher sensitivity in only 90 minutes with Human VEGF R1 ELISA Kit (FLT1) (**ab195210**) from our SimpleStep ELISA® range.

平台 Microplate

性能

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

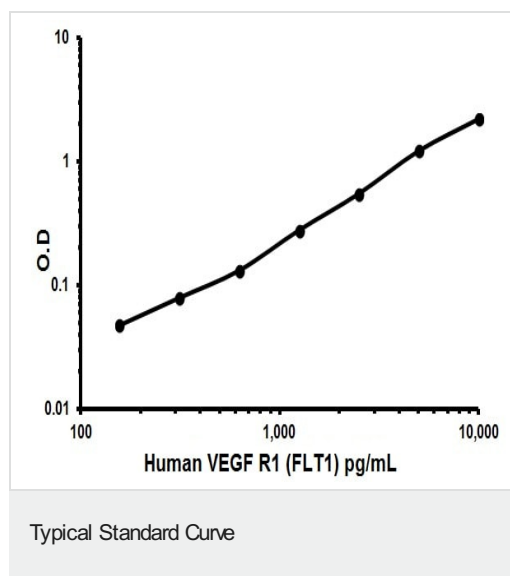
组件	标识符	1 x 96 tests
ABC Diluent Buffer	Blue Cap	1 x 12ml
Antibody Diluent Buffer	Green Cap	1 x 12ml
Anti-Human VEGF Receptor 1 antibody Microplate (12 x 8 wells)		1 unit
Avidin-Biotin-Peroxidase Complex (ABC)		1 x 100µl
Biotinylated anti-Human VEGF Receptor 1 antibody		1 x 100µl
Lyophilized recombinant Human VEGF Receptor 1 standard		2 x 10ng
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 Receptor for VEGF, VEGFB and PGF. Has a tyrosine-protein kinase activity. The VEGF-kinase ligand/receptor signaling system plays a key role in vascular development and regulation of vascular permeability. Isoform sFlt1 may have an inhibitory role in angiogenesis.

组织特异性 Mostly in normal lung, but also in placenta, liver, kidney, heart and brain tissues. Specifically expressed in most of the vascular endothelial cells, and also expressed in peripheral blood monocytes. Isoform sFlt1 is strongly expressed in placenta.

序列相似性 Belongs to the protein kinase superfamily. Tyr protein kinase family. CSF-1/PDGF receptor subfamily.
Contains 7 Ig-like C2-type (immunoglobulin-like) domains.
Contains 1 protein kinase domain.

细胞定位 Secreted and Cell membrane.



Representative standard curve using ab119613

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