

Human CD31 ELISA Kit (PECAM1) ab45910

1 References 1 图像

概述

产品名称	人CD31 ELISA试剂盒(PECAM1)				
检测方法	Colorimetric				
精确度	批次内				
	样品	n	Mean	SD	CV%
	Sample A	11	5.323ng/ml	0.19	3.5%
	Sample B	11	1.236ng/ml	0.077	6.2%
	批次间				
	样品	n	Mean	SD	CV%
	Sample A	11	5.406ng/ml	0.292	5.4%
	Sample B	11	1.205ng/ml	0.058	4.8%
样品类型	Cell culture supernatant, Serum, Plasma				
检测类型	Sandwich (quantitative)				
灵敏度	< 0.04 ng/ml				
范围	0.31 ng/ml - 10 ng/ml				
检测时间	2h 45m				
实验步骤	Multiple steps standard assay				
种属反应性	与反应: Human				
产品概述	Abcam's CD31 (PECAM1) Human <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of CD31 in Human serum, plasma, buffered solutions and cell culture medium				

A monoclonal antibody specific for CD31 has been coated onto the wells of the microtiter strips provided. Samples, including standards of known CD31 concentrations, control specimens or unknowns are pipetted into these wells. During the first incubation, the standards or samples and a biotinylated monoclonal antibody specific for CD31 are simultaneously incubated. After washing, the enzyme Streptavidin-HRP, that binds the biotinylated antibody is added, incubated

and washed. A TMB substrate solution is added which acts on the bound enzyme to induce a colored reaction product. The intensity of this colored product is directly proportional to the concentration of CD31 present in the samples.

This kit will recognize both endogenous and recombinant Human CD31.

Get results in 90 minutes with Human CD31 ELISA Kit (PECAM) ([ab190814](#)) from our SimpleStep ELISA® range.

平台

Microplate

性能

存放说明

Store at +4°C. Please refer to protocols.

组件	标识符	1 x 96 tests	2 x 96 tests
10X Standard Diluent Buffer	Black	1 x 15ml	1 x 25ml
200X Wash Buffer	White	1 x 10ml	2 x 10ml
Biotinylated Antibody Diluent	Red	1 x 7ml	1 x 13ml
Biotinylated anti-CD31 antibody	Red	1 x 400µl	2 x 400µl
CD31 Control	Silver	2 vials	4 vials
CD31 Microplate (12 x 8 well strips)		1 unit	2 units
CD31 Standard	Yellow	2 vials	4 vials
Chromogen TMB Substrate Solution		1 x 11ml	1 x 24ml
HRP Diluent	Red	1 x 12ml	1 x 23ml
Stop Reagent	Black	1 x 11ml	2 x 11ml
Streptavidin-HRP		2 x 5µl	4 x 5µl

功能

Induces susceptibility to atherosclerosis (By similarity). Cell adhesion molecule which is required for leukocyte transendothelial migration (TEM) under most inflammatory conditions. Tyr-690 plays a critical role in TEM and is required for efficient trafficking of PECAM1 to and from the lateral border recycling compartment (LBRC) and is also essential for the LBRC membrane to be targeted around migrating leukocytes. Prevents phagocyte ingestion of closely apposed viable cells by transmitting 'detachment' signals, and changes function on apoptosis, promoting tethering of dying cells to phagocytes (the encounter of a viable cell with a phagocyte via the homophilic interaction of PECAM1 on both cell surfaces leads to the viable cell's active repulsion from the phagocyte. During apoptosis, the inside-out signaling of PECAM1 is somehow disabled so that the apoptotic cell does not actively reject the phagocyte anymore. The lack of this repulsion signal together with the interaction of the eat-me signals and their respective receptors causes the attachment of the apoptotic cell to the phagocyte, thus triggering the process of engulfment). Isoform Delta15 is unable to protect against apoptosis. Modulates BDKRB2 activation. Regulates

bradykinin- and hyperosmotic shock-induced ERK1/2 activation in human umbilical cord vein cells (HUVEC).

组织特异性

Expressed on platelets and leukocytes and is primarily concentrated at the borders between endothelial cells. Isoform Long predominates in all tissues examined. Isoform Delta12 is detected only in trachea. Isoform Delta14-15 is only detected in lung. Isoform Delta14 is detected in all tissues examined with the strongest expression in heart. Isoform Delta15 is expressed in brain, testis, ovary, cell surface of platelets, human umbilical vein endothelial cells (HUVECs), Jurkat T-cell leukemia, human erythroleukemia (HEL) and U937 histiocytic lymphoma cell lines (at protein level).

序列相似性

Contains 6 Ig-like C2-type (immunoglobulin-like) domains.

结构域

The Ig-like C2-type domains 2 and 3 contribute to formation of the complex with BDKRB2 and in regulation of its activity.

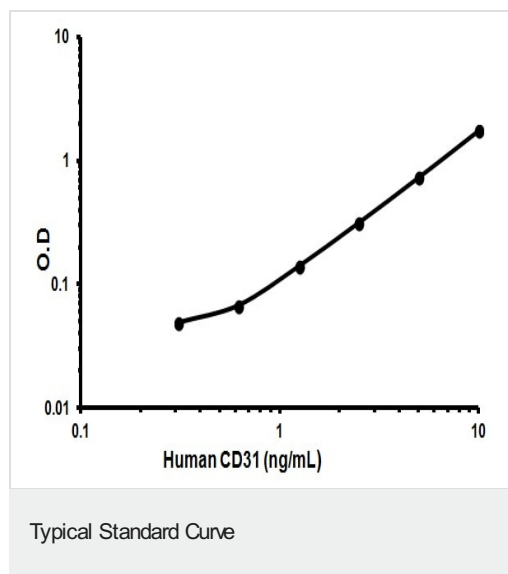
翻译后修饰

Phosphorylated on Ser and Tyr residues after cellular activation. In endothelial cells Fyn mediates mechanical-force (stretch or pull) induced tyrosine phosphorylation.

细胞定位

Membrane. Cell junction. Localizes to the lateral border recycling compartment (LBRC) and recycles from the LBRC to the junction in resting endothelial cells and Cell junction. Localizes to the lateral border recycling compartment (LBRC) and recycles from the LBRC to the junction in resting endothelial cells.

图片



Representative Standard Curve using ab45910.

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