



Product datasheet

anti-Bevacizumab ELISA Kit ab237658

1 图像

概述

产品名称	anti-Bevacizumab ELISA试剂盒				
检测方法	Colorimetric				
精确度	批次内				
	样品	n	Mean	SD	CV%
	Overall				< 15%
	批次间				
	样品	n	Mean	SD	CV%
	Overall				< 15%
样品类型	Serum, Plasma				
检测类型	Sandwich (quantitative)				
灵敏度	30 ng/ml				
范围	62 ng/ml - 500 ng/ml				
回收率	特定样本回收率				
	样品类型	平均%		范围	
	Serum			85% - 115%	
实验步骤	Multiple steps standard assay				
种属反应性	与反应: Human				
产品概述	anti-Bevacizumab ELISA Kit (ab237658) is designed to quantify/measure the antibody against Bevacizumab with high specificity and sensitivity in biological matrices.				
说明	This product is manufactured by BioVision, an Abcam company and was previously called E4389 BioSim™ anti-Bevacizumab (Human) ELISA Kit. E4389-100 is the same size as the 96 test size of ab237658.				
平台	Microplate				

性能

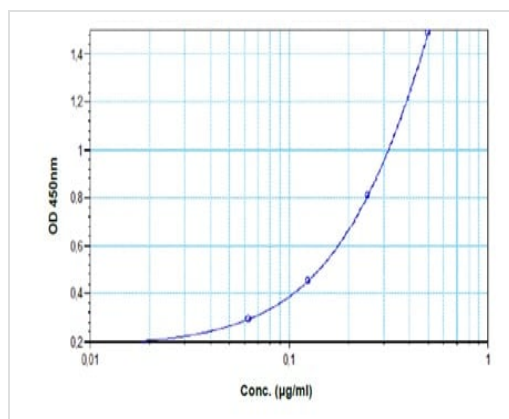
存放说明 Store at +4°C. Please refer to protocols.

组件	96 tests
Micro ELISA Plate	1 unit
Bevacizumab standard 1	1 x 1ml
Bevacizumab standard 2	1 x 1ml
Bevacizumab Standard 3	1 x 1ml
Bevacizumab Standard 4	1 x 1ml
Bevacizumab Standard 5	1 x 1ml
Bevacizumab Standard 6	1 x 1ml
Bevacizumab Standard 7	1 x 1ml
Assay Buffer	1 x 50ml
Confirmation Reagent	1 x 12ml
Peroxidase Conjugate	1 x 12ml
TMB substrate	1 x 12ml
Stop Solution	1 x 12ml
Wash buffer (20X)	1 x 50ml
Plate sealers	2 tablets

相关性

Bevacizumab is a recombinant human IgG1 monoclonal antibody specific for all human vascular endothelial growth factor-A (VEGF-A) isoforms. The humanized anti-VEGF monoclonal antibody, bevacizumab, has been approved by the FDA as a first-line treatment for metastatic colorectal cancer in combination with chemotherapy. The pharmacokinetic properties of bevacizumab in several species have been previously described and are consistent with a typical humanized monoclonal antibody. It was shown in the literature that the surveillance of circulating concentration during maintenance therapy represents a direct and/or indirect factor for some other side effects. Identification of biomarkers for (non-)response and risk factors for adverse drug reactions that might be related to serum concentrations and maintaining the effective concentration of Bevacizumab in order to potentially avoid some side effects with a reliable method might be beneficial.

图片



Typical standard curve.

Typical Standard Curve: This standard curve is for demonstration only. A standard curve must be run with each assay.

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