

Human MICA ELISA Kit ab59569

3 References 1 图像

概述

产品名称	人MICA ELISA试剂盒				
检测方法	Colorimetric				
精确度	批次内				
	样品	n	Mean	SD	CV%
	A	6	2214.6pg/ml	50.3	2.3%
	B	6	1102.1pg/ml	39.5	3.6%
	C	6	641.6pg/ml	47.7	7.4%
	批次间				
	样品	n	Mean	SD	CV%
	A	18	2416pg/ml	187.1	7.7%
	B	16	1200pg/ml	105.2	8.8%
	C	18	627pg/ml	50.9	8.1%
样品类型	Cell culture supernatant				
检测类型	Sandwich (quantitative)				
灵敏度	= 123 pg/ml				
范围	156.25 pg/ml - 5000 pg/ml				
检测时间	3h 45m				
实验步骤	Multiple steps standard assay				
种属反应性	与反应: Human				
产品概述	Abcam's MICA Human <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for				

unknowns are pipetted into these wells. During the first incubation, the standards or samples and a biotinylated monoclonal antibody specific for MICA 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 MICA present in the samples.

This kit will recognize both endogenous and recombinant Human MICA.

平台 Microplate

性能

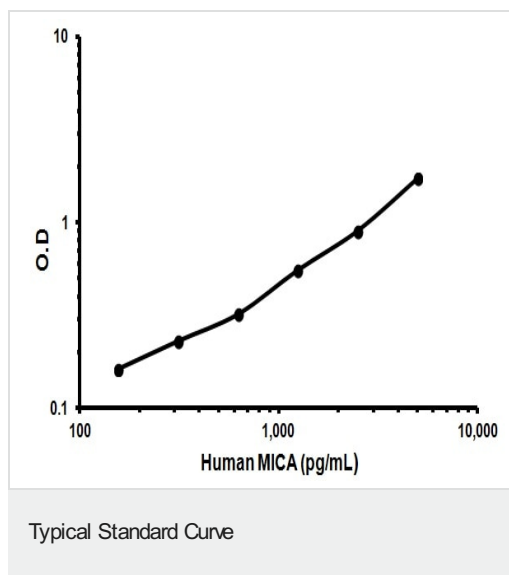
存放说明 Store at +4°C. Please refer to protocols.

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

相关性 The MHC class I chain-related (MIC) proteins are related to the Major histocompatibility complex (MHC) class I proteins which are ubiquitously expressed and mediate the recognition of intracellular antigens by cytotoxic T cells. The MHC class I chain-related (MIC) proteins are recognized by NKG2D, a receptor on NK and T cells, and promote anti-tumor activity. MICA, a member of the MIC family, is widely expressed on many tumors, and it is the MICA/NKG2D interaction that is thought to stimulate the anti-tumor reactivity by T lymphocytes. MICA is present in virtually every tissue except the nervous system, suggesting that MIC protein expression may only be one component of the anti-tumor activity of the immune system.

细胞定位 Plasma membrane

图片



Representative Standard Curve using ab59569.

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