

# Human MICA ELISA Kit ab100592

2 References 3 图像

概述

产品名称	人MICA ELISA试剂盒
检测方法	Colorimetric
样品类型	Cell culture supernatant, Serum, Plasma
检测类型	Sandwich (quantitative)
灵敏度	< 20 pg/ml
范围	13.72 pg/ml - 10000 pg/ml
回收率	100 %

特定样本回收率

样品类型	平均 %	范围
Cell culture supernatant	116.9	109% - 124%
Serum	112.3	95% - 129%
Plasma	93.83	75% - 107%

实验步骤	Multiple steps standard assay
种属反应性	与反应: Human
产品概述	<p>Abcam's MICA Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an <i>in vitro</i> enzyme linked immunosorbent assay for the quantitative measurement of Human MICA in serum, plasma, and cell culture supernatants.</p> <p>This assay employs an antibody specific for Human MICA coated on a 96-well plate. Standards and samples are pipetted into the wells and MICA present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human MICA antibody is added. After washing away unbound biotinylated antibody, HRP-conjugated streptavidin is pipetted to the wells. The wells are again washed, a TMB substrate solution is added to the wells and color develops in proportion to the amount of MICA bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.</p>
说明	Optimization may be required with urine samples.
平台	Microplate

性能

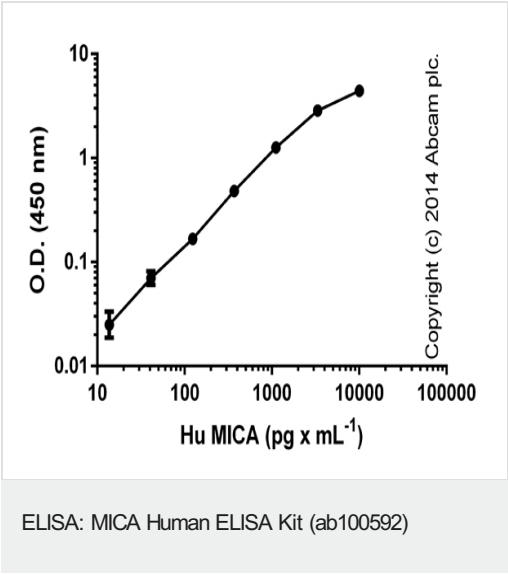
存放说明 Store at -20°C. Please refer to protocols.

组件	1 x 96 tests
20X Wash Buffer	1 x 25ml
300X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent	1 x 15ml
Biotinylated anti-Human MICA	2 vials
MICA Microplate (12 x 8 wells)	1 unit
Recombinant Human MICA Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

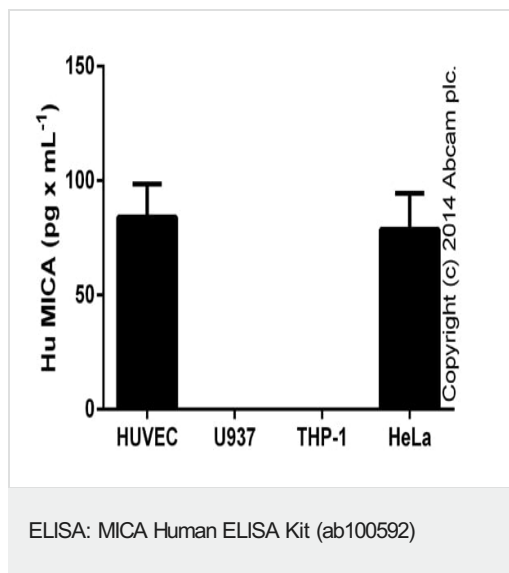
**相关性** The MHC class I chain-related (MIC) proteins are related to the Major histocompatibility complex (MHC) class I proteins which are ubiquitously expressed and mediate the recognition of intracellular antigens by cytotoxic T cells. The MHC class I chain-related (MIC) proteins are recognized by NKG2D, 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

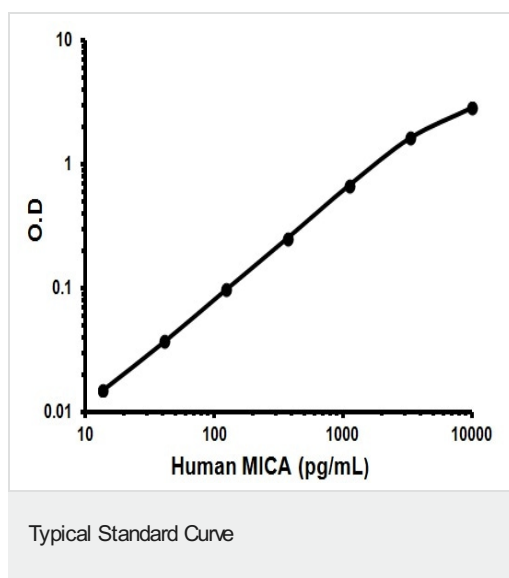
图片



Standard curve with background signal subtracted (duplicates; +/- SD).



MICA measured in undiluted cell culture supernatants, U937 and THP-1 signals were below level of detection (13.7 pg x mL<sup>-1</sup>) (duplicates +/- SD).



Representative Standard Curve using ab100592.

**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