

Human MICB ELISA Kit ab100593

2 References 画像数 2

製品の概要

製品名	Human MICB ELISA Kit
検出方法	Colorimetric
サンプルの種類	Cell culture supernatant, Serum, Plasma
アッセイタイプ	Sandwich (quantitative)
検出感度	< 75 pg/ml
検出範囲	0.069 ng/ml - 50 ng/ml
添加回収試験	95 %

特定サンプルでの回収試験

サンプルの種類	平均 %	測定範囲
Cell culture supernatant	98.22	91% - 125%
Serum	113.1	102% - 122%
Plasma	75.47	65% - 88%

ステップ Multiple steps standard assay

種交差性 交差種: Human

製品の概要 Abcam's MICB Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme linked immunosorbent assay for the quantitative measurement of Human MICB in serum, plasma, and cell culture supernatants.

This assay employs an antibody specific for Human MICB coated on a 96-well plate. Standards and samples are pipetted into the wells and MICB present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human MICB 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 MICB bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

特記事項 Optimization may be required with urine samples.

試験プラットフォーム Microplate

## 製品の特性

### 保存方法

Store at -20°C. Please refer to protocols.

内容	1 x 96 tests
200X HRP-Streptavidin Concentrate	1 x 200µl
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
Assay Diluent A	1 x 30ml
Biotinylated anti-Human MICB	2 vials
MICB Microplate (12 x 8 wells)	1 unit
Recombinant Human MICB Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

### 機能

Seems to have no role in antigen presentation. Acts as a stress-induced self-antigen that is recognized by gamma delta T cells. Ligand for the KLRK1/NKG2D receptor. Binding to KLRK1 leads to cell lysis.

### 組織特異性

Widely expressed with the exception of the central nervous system where it is absent. Expressed in many, but not all, epithelial tumors of lung, breast, kidney, ovary, prostate and colon. In hepatocellular carcinomas, expressed in tumor cells but not in surrounding non-cancerous tissue.

### 関連疾患

Genetic variations in MICA are a cause of susceptibility to rheumatoid arthritis (RA) [MIM:180300]. It is a systemic inflammatory disease with autoimmune features and a complex genetic component. It primarily affects the joints and is characterized by inflammatory changes in the synovial membranes and articular structures, widespread fibrinoid degeneration of the collagen fibers in mesenchymal tissues, and by atrophy and rarefaction of bony structures.  
Note=The MICB\*004 allele is associated with rheumatoid arthritis.  
Note=Genetic variation in MICB is associated with cytomegalovirus and herpes simplex virus I seropositivity and this may be associated with schizophrenia risk.

### 配列類似性

Belongs to the MHC class I family. MIC subfamily.  
Contains 1 Ig-like C1-type (immunoglobulin-like) domain.

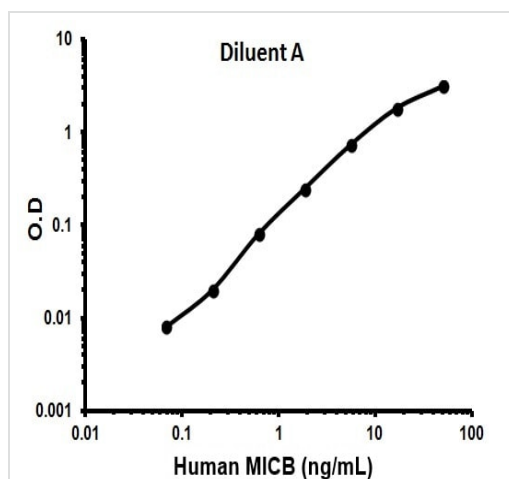
### 翻訳後修飾

Proteolytically cleaved and released from the cell surface of tumor cells.

### 細胞内局在

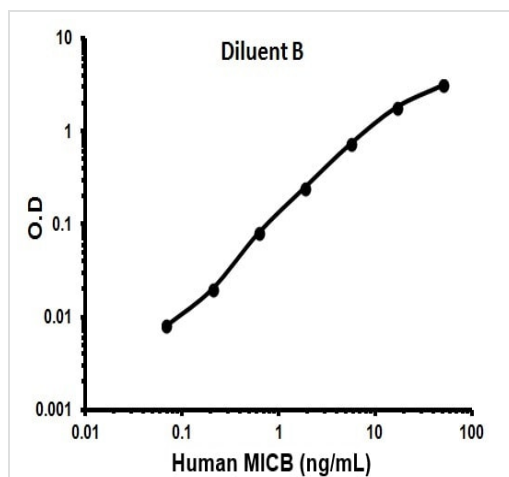
Cell membrane. Binding to human cytomegalovirus glycoprotein UL16 causes sequestration in the endoplasmic reticulum.

## 画像



Representative Standard Curve using ab100593.

Typical Standard Curve



Representative Standard Curve using ab100593.

Typical Standard Curve

**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.co.jp/abpromise> or contact our technical team.

## Terms and conditions

---

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors