

Human MICB ELISA Kit ab267804

重组 SimpleStep ELISA

4 图像

概述

产品名称 人MICB ELISA试剂盒

检测方法 Colorimetric

精确度

批次内

样品	n	Mean	SD	CV%
Serum	8			3.8%

批次间

样品	n	Mean	SD	CV%
Serum	3			2.5%

样品类型 Cell culture supernatant, Serum, Hep Plasma, EDTA Plasma, Cit plasma

检测类型 Sandwich (quantitative)

灵敏度 59.9 pg/ml

范围 156.25 pg/ml - 10000 pg/ml

回收率

特定样本回收率

样品类型	平均%	范围
Cell culture supernatant	115	112% - 117%
Serum	100	100% - 100%
Hep Plasma	90	89% - 90%
EDTA Plasma	97	97% - 98%
Cit plasma	97	96% - 98%

检测时间 1h 30m

实验步骤 One step assay

种属反应性

与反应: Human

产品概述

Human MICB ELISA Kit (ab267804) is a single-wash 90 min sandwich ELISA designed for the quantitative measurement of MICB protein in cell culture supernatant, edta plasma, hep plasma, serum, and cit plasma. It uses our proprietary SimpleStep ELISA® technology. Quantitate Human MICB with 59.9 pg/ml sensitivity.

SimpleStep ELISA® technology employs capture antibodies conjugated to an affinity tag that is recognized by the monoclonal antibody used to coat our SimpleStep ELISA® plates. This approach to sandwich ELISA allows the formation of the antibody-analyte sandwich complex in a single step, significantly reducing assay time. See the SimpleStep ELISA® protocol summary in the image section for further details. Our SimpleStep ELISA® technology provides several benefits:

- Single-wash protocol reduces assay time to 90 minutes or less
- High sensitivity, specificity and reproducibility from superior antibodies
- Fully validated in biological samples
- 96-wells plate breakable into 12 x 8 wells strips

A 384-well SimpleStep ELISA® microplate (**ab203359**) is available to use as an alternative to the 96-well microplate provided with SimpleStep ELISA® kits.

说明

MHC class I polypeptide-related sequence B (MICB) is a stress induced ligand for the KLRK1/NKG2D receptor that neither associate with β_2 -microglobins nor bind peptides. The MICB protein is encoded by the *MICB* gene located within the MHC locus and are commonly expressed on tumors of epithelial origin.

平台

Pre-coated microplate (12 x 8 well strips)

性能**存放说明**

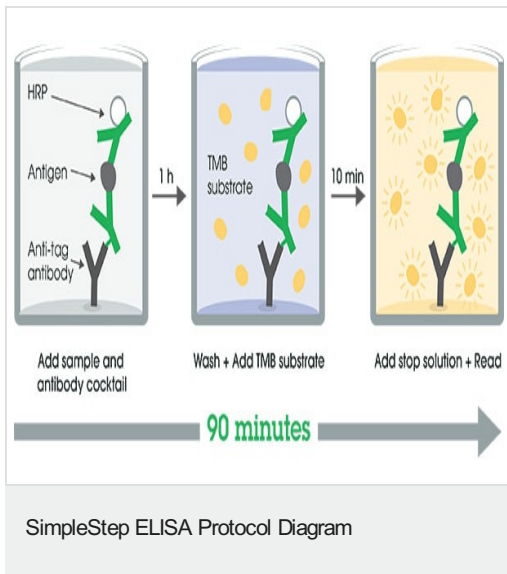
Store at +4°C. Please refer to protocols.

组件	1 x 96 tests
10X Human MICB Capture Antibody	1 x 600µl
10X Human MICB Detector Antibody	1 x 600µl
10X Wash Buffer PT (ab206977)	1 x 20ml
Antibody Diluent CPI2	1 x 6ml
Human MICB Lyophilized Recombinant Protein	2 vials
Plate Seals	1 unit
Sample Diluent NS (ab193972)	1 x 50ml
SimpleStep Pre-Coated 96-Well Microplate (ab206978)	1 unit

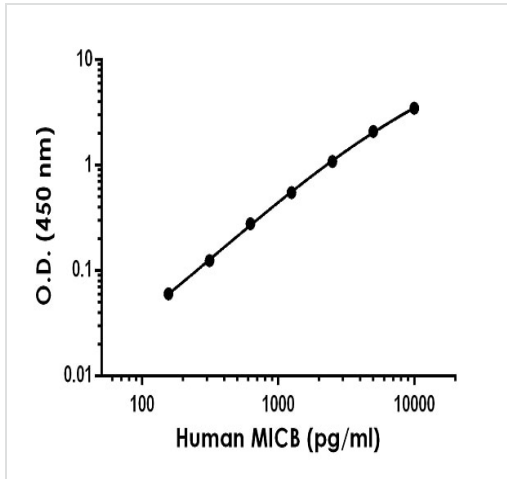
组件	1 x 96 tests
Stop Solution	1 x 12ml
TMB Development Solution	1 x 12ml

功能	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.
组织特异性	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.
疾病相关	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.
序列相似性	Belongs to the MHC class I family. MIC subfamily. Contains 1 Ig-like C1-type (immunoglobulin-like) domain.
翻译后修饰	Proteolytically cleaved and released from the cell surface of tumor cells.
细胞定位	Cell membrane. Binding to human cytomegalovirus glycoprotein UL16 causes sequestration in the endoplasmic reticulum.

图片

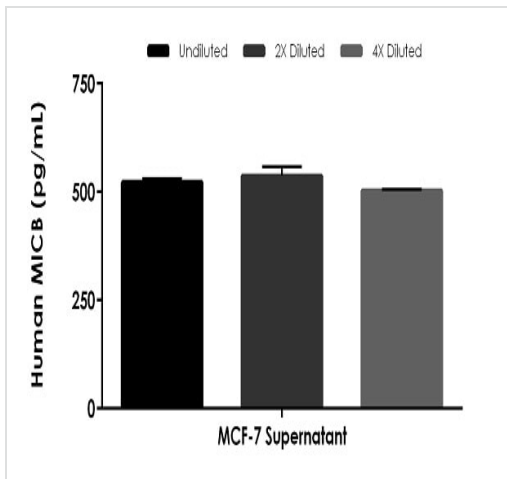


SimpleStep ELISA technology allows the formation of the antibody-antigen complex in one single step, reducing assay time to 90 minutes. Add samples or standards and antibody mix to wells all at once, incubate, wash, and add your final substrate. See protocol for a detailed step-by-step guide.



Example of human MICB standard curve in Sample Diluent NS.

The MICB standard curve was prepared as described in Section 10. Raw data values are shown in the table. Background-subtracted data values (mean +/- SD) are graphed.



Interpolated concentrations of native MICB in human cell culture supernatant sample.

The concentrations of MICB were measured in duplicates, interpolated from the MICB standard curves and corrected for sample dilution. Undiluted sample is as follows: MCF-7 supernatant 100%. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean MICB concentration was determined to be 520.4 pg/mL in MCF-7 supernatant.

Powered by
recombinant antibodies



Research with confidence
Consistent and reproducible results



Long-term and scalable supply
Recombinant technology



Success from the first experiment
Confirmed specificity



Ethical standards compliant
Animal-free production

Sandwich ELISA - Human MICB ELISA Kit
(ab267804)

To learn more about the advantages of recombinant antibodies see [here](#).

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