

Mouse KIM1 ELISA Kit (TIM-1) ab119596

10 References 1 图像

概述

产品名称

检测方法

精确度

小鼠KIM1 ELISA试剂盒(TIM-1)

Colorimetric

批次内

样品	n	Mean	SD	CV%
1	16	87pg/ml	6.87	= 7.9%
2	16	279pg/ml	14.78	= 5.3%
3	16	842pg/ml	36.2	= 4.3%

批次间

样品	n	Mean	SD	CV%
1	24	85pg/ml	7.31	= 8.6%
2	24	261pg/ml	14.87	= 5.7%
3	24	792pg/ml	38.8	= 4.9%

样品类型

检测类型

灵敏度

范围

实验步骤

种属反应性

产品概述

Cell culture supernatant, Urine, Serum, Hep Plasma, EDTA Plasma

Sandwich (quantitative)

< 2 pg/ml

31.2 pg/ml - 2000 pg/ml

Multiple steps standard assay

与反应: Mouse

Abcam's mouse KIM-1 (TIM-1) *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is

A KIM-1 (TIM-1) specific goat polyclonal antibody has been precoated onto 96-well plates. Standards and test samples are added to the wells and are incubated. A biotinylated detection

polyclonal antibody from goat specific for KIM-1 (TIM-1) is then added followed by washing with PBS or TBS buffer. Avidin-Biotin-Peroxidase Complex is added and unbound conjugates are washed away with PBS or TBS buffer. TMB is then used to visualize the HRP enzymatic reaction. TMB is catalyzed by HRP to produce a blue color product that changes into yellow after adding acidic stop solution. The density of yellow coloration is directly proportional to the mouse KIM-1 (TIM-1) amount of sample captured in plate.

Get higher sensitivity in only 90 minutes with Mouse KIM-1 ELISA Kit (TIM1) ([ab213477](#)) from our SimpleStep ELISA® range.

平台 Microplate

性能

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

组件	标识符	1 x 96 tests	1 x 96 tests
ABC Diluent Buffer	Blue Cap	1 x 12ml	1 x 12ml
Antibody Diluent Buffer	Green Cap	1 x 12ml	1 x 12ml
Anti-mouse KIM1 antibody Microplate (12 x 8 wells)		1 unit	1 unit
Avidin-Biotin-Peroxidase Complex (ABC)		1 x 100µl	1 x 100µl
Biotinylated anti-Mouse KIM1 antibody		1 x 100µl	1 x 100µl
Lyophilized recombinant Mouse KIM1 standard		2 x 10ng	2 x 10ng
Plate Seal		1 x 4 units	1 x 4 units
Sample Diluent Buffer	Green Cap	1 x 30ml	1 x 30ml
TMB Color Developing Agent	Amber Bottle	1 x 10ml	1 x 10ml
TMB Stop Solution	Yellow Cap	1 x 10ml	1 x 10ml
Wash Buffer (25X)		1 x 20ml	1 x 20ml

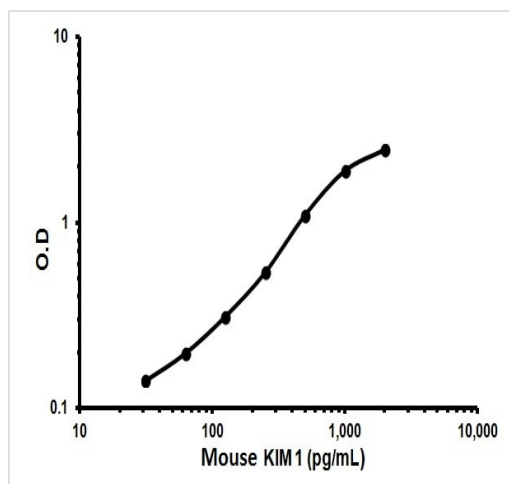
功能 May play a role in T-helper cell development and the regulation of asthma and allergic diseases. Receptor for TIMD4 (By similarity). In case of human hepatitis A virus (HHAV) infection, functions as a cell-surface receptor for the virus. May play a role in kidney injury and repair.

组织特异性 Widely expressed, with highest levels in kidney and testis. Expressed by activated CD4+ T-cells during the development of helper T-cells responses.

序列相似性 Belongs to the immunoglobulin superfamily. TIM family.
Contains 1 Ig-like V-type (immunoglobulin-like) domain.

细胞定位 Membrane.

图片



Representative Standard Curve for ab119596

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.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