

Mouse RAGE ELISA Kit ab100738

4 图像

概述

产品名称	小鼠RAGE 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	< 15 pg/ml
范围	13.72 pg/ml - 10000 pg/ml
回收率	87 %

特定样本回收率

样品类型	平均%	范围
Cell culture supernatant	75.89	67% - 82%
Serum	102.9	91% - 110%
Plasma	82.69	71% - 92%

实验步骤 Multiple steps standard assay

种属反应性 与反应: Mouse

产品概述 Abcam's RAGE Mouse ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of mouse RAGE in serum, plasma and cell culture supernatants.

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

平台 Microplate

性能

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

组件	1 x 96 tests
160X HRP-Streptavidin Concentrate	1 x 200µl
20X Wash Buffer	1 x 25ml
5X Assay Diluent	1 x 15ml
Biotinylated anti-Mouse RAGE	2 vials
RAGE Microplate (12 x 8 wells)	1 unit
Recombinant Mouse RAGE Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

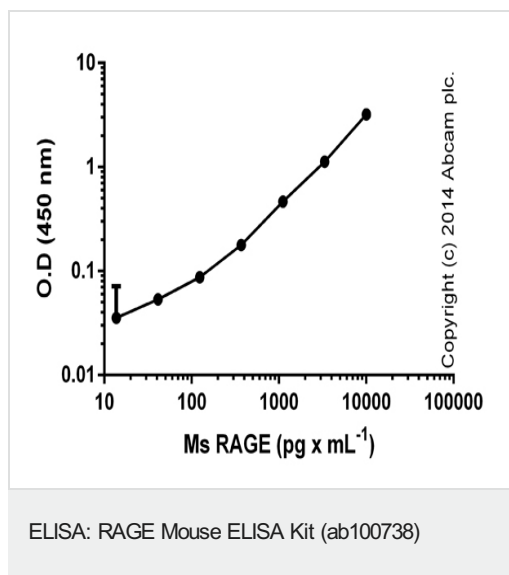
功能Mediates interactions of advanced glycosylation end products (AGE). These are nonenzymatically glycosylated proteins which accumulate in vascular tissue in aging and at an accelerated rate in diabetes. Acts as a mediator of both acute and chronic vascular inflammation in conditions such as atherosclerosis and in particular as a complication of diabetes. AGE/RAGE signaling plays an important role in regulating the production/expression of TNF-alpha, oxidative stress, and endothelial dysfunction in type 2 diabetes. Interaction with S100A12 on endothelium, mononuclear phagocytes, and lymphocytes triggers cellular activation, with generation of key proinflammatory mediators. Interaction with S100B after myocardial infarction may play a role in myocyte apoptosis by activating ERK1/2 and p53/TP53 signaling (By similarity). Receptor for amyloid beta peptide. Contributes to the translocation of amyloid-beta peptide (ABPP) across the cell membrane from the extracellular to the intracellular space in cortical neurons. ABPP-initiated RAGE signaling, especially stimulation of p38 mitogen-activated protein kinase (MAPK), has the capacity to drive a transport system delivering ABPP as a complex with RAGE to the intraneuronal space.

组织特异性Endothelial cells.

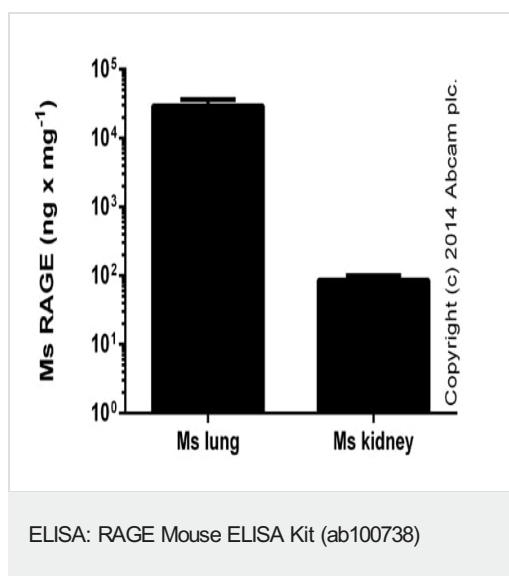
序列相似性Contains 2 Ig-like C2-type (immunoglobulin-like) domains.
Contains 1 Ig-like V-type (immunoglobulin-like) domain.

细胞定位Secreted and Cell membrane.

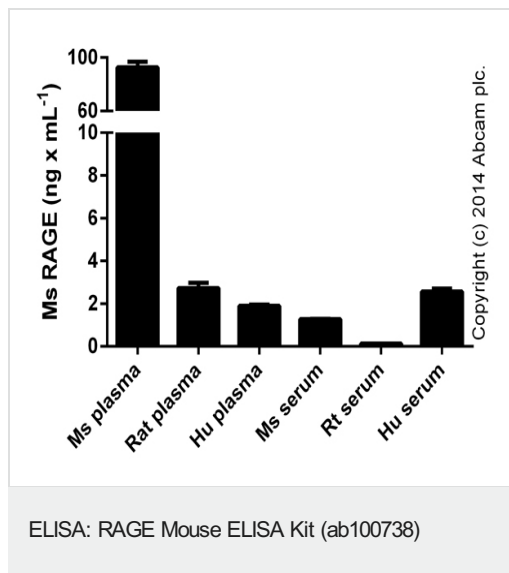
图片



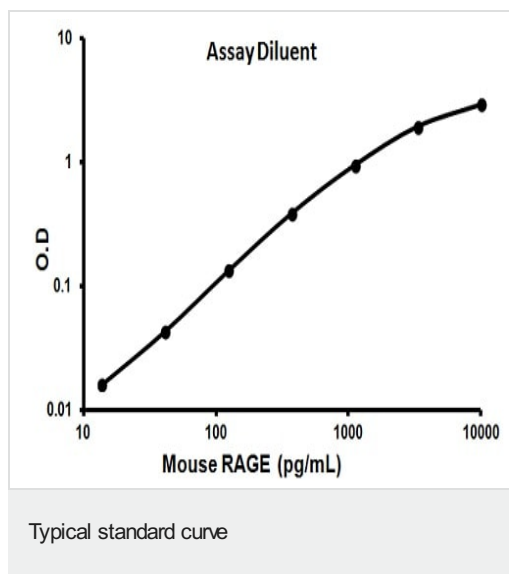
Standard curve of mouse RAGE with background signal subtracted (duplicates; \pm SD).



Mouse RAGE measured in mouse tissue lysates (in RIPA buffer tested in the range of 0.1 μg to 0.1 mg protein, quantity of RAGE shown as per mg of extracted protein), duplicates \pm SD.



Mouse RAGE measured in biological fluids (dilution range neat - 1/10; duplicates +/- SD).



Representative standard curve using ab100738

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