

Rat ICAM1 ELISA Kit (CD54) ab100763

9 References 2 图像

概述

| | |
|------|---|
| 产品名称 | 大鼠ICAM1 ELISA试剂盒(CD54) |
| 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 | 28.81 pg/ml - 7000 pg/ml |
| 回收率 | 91 % |

特定样本回收率

| 样品类型 | 平均% | 范围 |
|--------------------------|-------|------------|
| Cell culture supernatant | 98.54 | 89% - 107% |
| Serum | 89.42 | 80% - 101% |
| Plasma | 85.54 | 76% - 95% |

实验步骤 Multiple steps standard assay

种属反应性 与反应: Rat

产品概述 Abcam's ICAM1 Rat ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Rat ICAM- 1 in serum, plasma and cell culture supernatants.

This assay employs an antibody specific for Rat ICAM1 coated on a 96- well plate. Standards and samples are pipetted into the wells and ICAM1 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Rat ICAM1 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 ICAM1 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.

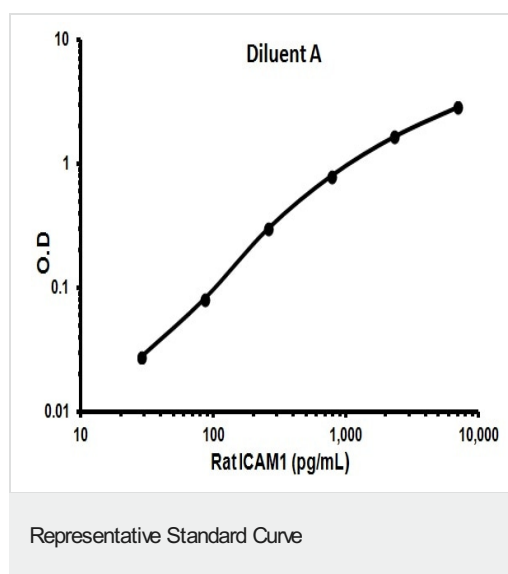
| 组件 | 1 x 96 tests |
|--|--------------|
| 20X Wash Buffer | 1 x 25ml |
| 400X HRP-Streptavidin Concentrate | 1 x 200µl |
| 5X Assay Diluent B | 1 x 15ml |
| Assay Diluent A | 1 x 30ml |
| Biotinylated anti-Rat ICAM1 | 2 vials |
| ICAM1 Microplate (12 x 8 wells) | 1 unit |
| Recombinant Rat ICAM1 Standard (lyophilized) | 2 vials |
| Stop Solution | 1 x 8ml |
| TMB One-Step Substrate Reagent | 1 x 12ml |

相关性

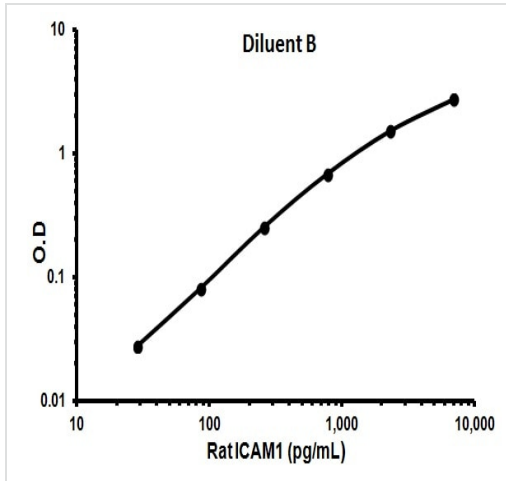
ICAM1 is a 85-110 kDa single chain type 1 integral membrane glycoprotein with an extracellular domain of five immunoglobulin superfamily repeats, a transmembrane region and a cytoplasmic domain. It shares considerable amino acid sequence homology with ICAM3 and with ICAM2. ICAM1 is expressed by activated endothelial cells. It is detected on cells of many other lineages (e.g. epithelial cells, fibroblasts, chondrocytes, B lymphocytes, T lymphocytes (low), monocytes, macrophages, dendritic cells and neutrophils), with lower levels that increase in inflammation. ICAM1 is also detected in some carcinoma and melanoma cells. Soluble ICAM1 is detectable in the plasma and is elevated in patients with various inflammatory syndromes. It is the receptor for rhinoviruses and malaria.

细胞定位

Membrane

图片

Representative Standard Curve using ab100763



Representative Standard Curve using ab100763

Representative 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