

Mouse ICAM1 ELISA Kit (CD54) ab100688

13 References 1 Image

Overview

|                  |   |
|------------------|---|
| Product name     | Mouse ICAM1 ELISA Kit (CD54)            |
| Detection method | Colorimetric                            |
| Sample type      | Cell culture supernatant, Serum, Plasma |
| Assay type       | Sandwich (quantitative)                 |
| Sensitivity      | < 20 pg/ml                              |
| Range            | 24.69 pg/ml - 6000 pg/ml                |
| Recovery         | 97 %                                    |

Sample specific recovery

| Sample type              | Average % | Range      |
|--------------------------|-----------|------------|
| Cell culture supernatant | 93.54     | 84% - 103% |
| Serum                    | 94.56     | 84% - 103% |
| Plasma                   | 115.2     | 97% - 127% |

|                    |   |
|--------------------|---|
| Assay duration     | Multiple steps standard assay   |
| Species reactivity | <b>Reacts with:</b> Mouse   |
| Product overview   | <p>Abcam's ICAM1 Mouse ELISA (Enzyme-Linked Immunosorbent Assay) kit is an <i>in vitro</i> enzyme-linked immunosorbent assay for the quantitative measurement of Mouse ICAM1 in serum, plasma and cell culture supernatants.</p> <p>This assay employs an antibody specific for mouse 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-mouse 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.</p> |
| Platform           | Microplate  |

Properties

**Storage instructions**

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

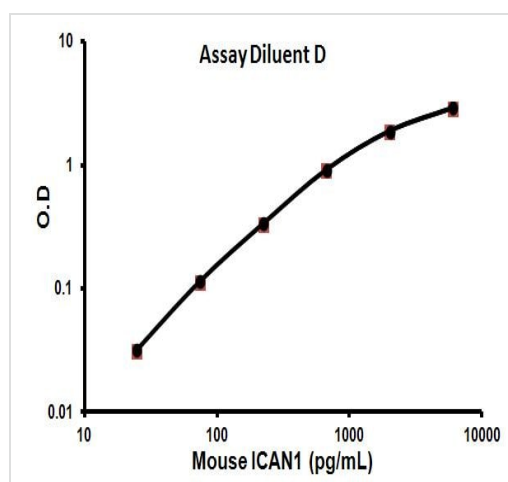
| Components                                     | 1 x 96 tests |
|--|--------------|
| 20X Wash Buffer                                | 1 x 25ml     |
| 400X HRP-Streptavidin Concentrate              | 1 x 200µl    |
| 5X Assay Diluent B                             | 1 x 15ml     |
| 5X Assay Diluent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

**Relevance**

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.

**Cellular localization**

Membrane

**Images**

Typical Standard Curve

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.com/abpromise> or contact our technical team.

### **Terms and conditions**

---

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