

Mouse ICAM1 ELISA Kit (CD54) ab100688

13 References 画像数 1

製品の概要

製品名	Mouse ICAM1 ELISA Kit (CD54)
検出方法	Colorimetric
サンプルの種類	Cell culture supernatant, Serum, Plasma
アッセイタイプ	Sandwich (quantitative)
検出感度	< 20 pg/ml
検出範囲	24.69 pg/ml - 6000 pg/ml
添加回収試験	97 %

特定サンプルでの回収試験

サンプルの種類	平均 %	測定範囲
Cell culture supernatant	93.54	84% - 103%
Serum	94.56	84% - 103%
Plasma	115.2	97% - 127%

ステップ	Multiple steps standard assay
種交差性	交差種: Mouse

製品の概要	<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> <p>試験プラットフォーム</p> <p>Microplate</p>
-------	---

製品の特性

保存方法

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

内容	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

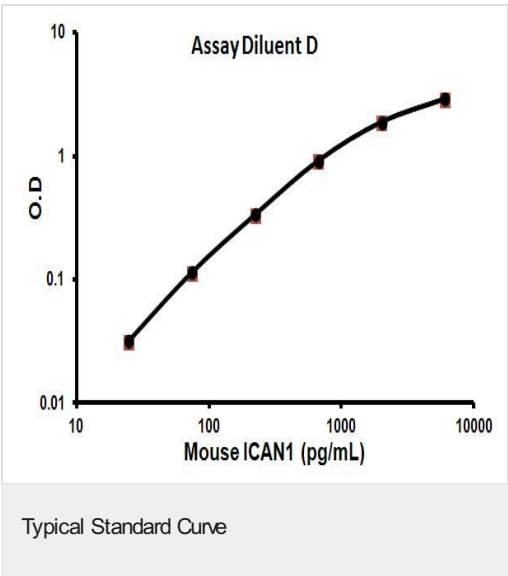
関連性

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.

細胞内局在

Membrane

画像



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.co.jp/abpromise> or contact our technical team.

Terms and conditions

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