

Human ICAM1 ELISA Set (without plates) ab47349

3 References 1 Image

Overview

Product name	Human ICAM1 ELISA Set (without plates)
Sample type	Cell culture supernatant, Serum, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	< 0.1 ng/ml
Range	0.25 ng/ml - 8 ng/ml
Assay time	1h 45m
Assay duration	Multiple steps standard assay
Species reactivity	<b>Reacts with:</b> Human
Product overview	<p>Abcam's ICAM1 Human <i>in vitro</i> ELISA Set is designed for the quantitative measurement of ICAM1 in cell culture supernatants, buffered solutions, serum, plasma and other body fluids.</p> <p>This ELISA set is intended to measure human ICAM1 in cell -culture supernatants, as well as in serum and plasma samples. The enclosed procedures provide general guidelines for ELISA preparation. This procedure may need some optimization depending on the plastic, matrix, standard range.</p> <p>The ICAM1 ELISA Set includes both the capture and the detection antibodies as well as Horse-Radish peroxidase conjugated Streptavidin (HRP-Strep) and TMB which are the reagents used in the colour development step.</p> <p>Recognizes both natural and recombinant human ICAM1.</p>
Tested applications	<b>Suitable for:</b> Sandwich ELISA
Platform	Reagents

Properties

Storage instructions Store at +4°C. Please refer to protocols.

Components	10 x 96 tests	5 x 96 tests	15 x 96 tests
Chromogen TMB Substrate Solution	4 vials	2 vials	6 vials

Components	10 x 96 tests	5 x 96 tests	15 x 96 tests
Detection biotinylated anti-ICAM1 antibody	2 x 0.55ml	1 x 0.55ml	3 x 0.55ml
ICAM1 Capture Antibody	2 x 0.5ml	1 x 0.5ml	3 x 0.5ml
ICAM1 Standard	10 vials	5 vials	15 vials
Streptavidin-HRP	2 vials	1 vial	3 vials

#### Function

ICAM proteins are ligands for the leukocyte adhesion protein LFA-1 (integrin alpha-L/beta-2). During leukocyte trans-endothelial migration, ICAM1 engagement promotes the assembly of endothelial apical cups through ARHGEF26/SGEF and RHOG activation. In case of rhinovirus infection acts as a cellular receptor for the virus.

#### Sequence similarities

Belongs to the immunoglobulin superfamily. ICAM family.  
Contains 5 Ig-like C2-type (immunoglobulin-like) domains.

#### Post-translational modifications

Monoubiquitinated, which is promoted by MARCH9 and leads to endocytosis.

#### Cellular localization

Membrane.

### Applications

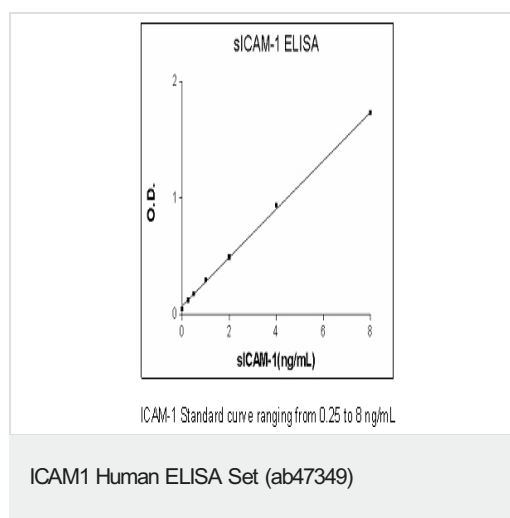
#### The Abpromise guarantee

Our [Abpromise guarantee](#) covers the use of ab47349 in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

Application	Abreviews	Notes
<b>Sandwich ELISA</b>		Use at an assay dependent concentration.

### Images



**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