

Human ICAM1 ELISA Set (without plates) ab47349

3 References 1 图像

概述

产品名称	人ICAM1 ELISA试剂盒(without plates)
样品类型	Cell culture supernatant, Serum, Plasma
检测类型	Sandwich (quantitative)
灵敏度	< 0.1 ng/ml
范围	0.25 ng/ml - 8 ng/ml
检测时间	1h 45m
实验步骤	Multiple steps standard assay
种属反应性	与反应: Human
产品概述	<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>
经测试应用	适用于: Sandwich ELISA
平台	Reagents

性能

存放说明 Store at +4°C. Please refer to protocols.

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

组件	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

**功能** 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 RHO G activation. In case of rhinovirus infection acts as a cellular receptor for the virus.

**序列相似性** Belongs to the immunoglobulin superfamily. ICAM family.  
Contains 5 Ig-like C2-type (immunoglobulin-like) domains.

**翻译后修饰** Monoubiquitinated, which is promoted by MARCH9 and leads to endocytosis.

**细胞定位** Membrane.

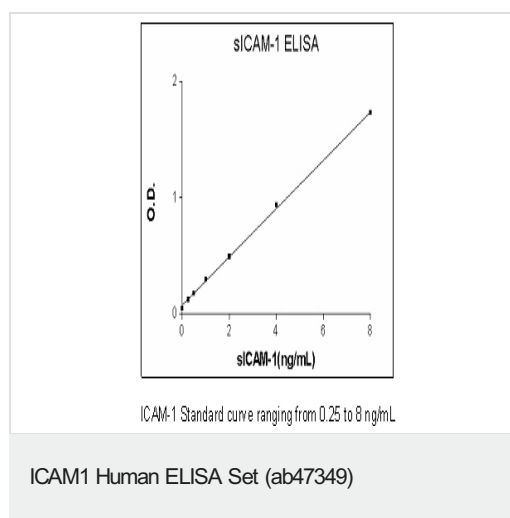
## 应用

**The Abpromise guarantee** **Abpromise™**承诺保证使用ab47349于以下的经测试应用

“应用说明”部分 下显示的仅为推荐的起始稀释度;实际最佳的稀释度/浓度应由使用者检定。

应用	Ab评论	说明
<b>Sandwich ELISA</b>		Use at an assay dependent concentration.

## 图片



**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