

Product datasheet

Human ICAM2 Matched Antibody Pair Kit ab217614

Recombinant

2 Images

Overview

<b>Product name</b>	Human ICAM2 Matched Antibody Pair Kit
<b>Detection method</b>	Colorimetric
<b>Assay type</b>	ELISA set
<b>Sensitivity</b>	9.17 pg/ml
<b>Range</b>	187.5 pg/ml - 12000 pg/ml
<b>Species reactivity</b>	<b>Reacts with:</b> Human
<b>Product overview</b>	<p>Human ICAM2 Matched Antibody Pair Kits include a capture and a biotinylated detector antibody pair, along with a calibrated protein standard, suitable for sandwich ELISA. The Matched Antibody Pair Kit can be used to quantify native and recombinant human ICAM2.</p> <p>Optimization of the kit reagents to sample type, immunoassay format or instrumentation may be required. Guidelines for use of this kit in a standard 96-well microplate sandwich ELISA using HRP/TMB system of colorimetric detection is described in this assay procedure for the purposes of quantification.</p> <p>Protocol information and tips on the use of the Matched Antibody Pair kits for sandwich ELISA can be found on our <a href="#">website</a>. An accessory pack can be purchased which includes buffer reagents required to perform 10 x 96-well plate sandwich ELISAs (<a href="#">ab210905</a>).</p> <p>For additional information on the performance of the antibody pair used in this kit, please see our equivalent SimpleStep ELISA kit <a href="#">ab212164</a>. Please note that while the antibody pair is the same provided in the corresponding SimpleStep ELISA Kit, due to differences in their formulation, this antibody pair cannot be used with the consumables provided with our SimpleStep ELISA Kits.</p>
<b>Tested applications</b>	<b>Suitable for:</b> ELISA
<b>Platform</b>	Reagents

Properties

**Storage instructions** Store at -20°C. Please refer to protocols.

Components	10 x 96 tests	5 x 96 tests
Human ICAM2 Capture Antibody	2 x 50µg	1 x 50µg
Human ICAM2 Detector Antibody	2 x 12.5µg	1 x 12.5µg
Human ICAM2 Lyophilized Protein	2 vials	1 vial

## Relevance

ICAM2 is a member of the intercellular adhesion molecule (ICAM) family. All ICAM proteins are type I transmembrane glycoproteins, contain 2-9 immunoglobulin like C2 type domains, and bind to the leukocyte adhesion LFA1 protein. This protein may play a role in lymphocyte recirculation by blocking LFA1 dependent cell adhesion. It mediates adhesive interactions important for antigen specific immune response, NK cell mediated clearance, lymphocyte recirculation, and other cellular interactions important for immune response and surveillance.

## Cellular localization

Membrane; Single-pass type I membrane protein

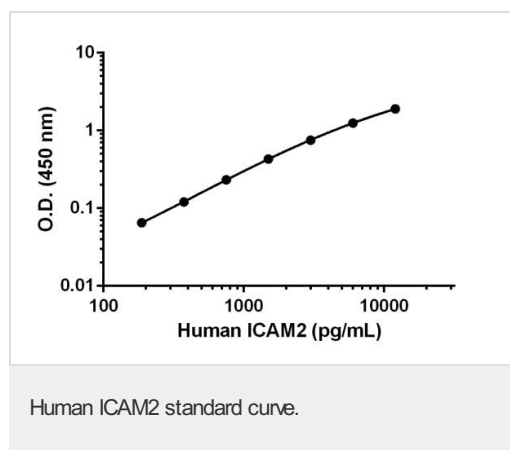
## Applications

Our [Abpromise guarantee](#) covers the use of **ab217614** 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
ELISA		Use at an assay dependent concentration.

## Images



Standard calibration curve. Background subtracted values are graphed.

Powered by  
recombinant antibodies



**Research with confidence**  
Consistent and reproducible results



**Long-term and scalable supply**  
Recombinant technology



**Success from the first experiment**  
Confirmed specificity



**Ethical standards compliant**  
Animal-free production

Sandwich ELISA - Human ICAM2 Matched Antibody  
Pair Kit (ab217614)

To learn more about the advantages of recombinant antibodies see [here](#).

**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