

Human ICAM3 ELISA Kit (CD50) ab46024

1 Image

Overview

Product name Human ICAM3 ELISA Kit (CD50)

Detection method Colorimetric

Precision	Intra-assay				
	Sample	n	Mean	SD	CV%
	Sample A	8	19.75ng/ml	0.172	0.86%
	Sample B	8	2.83ng/ml	0.099	3.49%

	Inter-assay				
	Sample	n	Mean	SD	CV%
	Sample A	36	19.51ng/ml	0.391	1.99%
	Sample B	36	2.71ng/ml	0.176	6.27%

Sample type Cell culture supernatant, Serum, Plasma

Assay type Sandwich (quantitative)

Sensitivity < 0.3 ng/ml

Range 0.62 ng/ml - 20 ng/ml

Assay time 1h 45m

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview

Abcam's ICAM3 (CD50) Human *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of ICAM3 in serum, plasma, buffered solutions and cell culture medium.

A monoclonal antibody specific for ICAM3 has been coated onto the wells of the microtiter strips provided. Samples, including standards of known ICAM3 concentrations, control specimens or unknowns are pipetted into these wells. During the first incubation, the standards or samples and a biotinylated monoclonal antibody specific for ICAM3 are simultaneously incubated. After washing, the enzyme Streptavidin-HRP, that binds the biotinylated antibody is added, incubated and washed. A TMB substrate solution is added which acts on the bound enzyme to induce a

colored reaction product. The intensity of this colored product is directly proportional to the concentration of ICAM3 present in the samples.

This kit will recognize both endogenous and recombinant Human ICAM3.

Platform

Microplate

Properties

Storage instructions

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

Components	Identifier	1 x 96 tests	2 x 96 tests
10X Standard Diluent Buffer	Black	1 x 15ml	1 x 25ml
200X Wash Buffer	White	1 x 10ml	2 x 10ml
Biotinylated Antibody Diluent	Red	1 x 7.5ml	1 x 13ml
Biotinylated anti-ICAM-3 (CD50)	Red	1 x 0.4ml	2 x 0.4ml
Chromogen TMB Substrate Solution		1 x 11ml	1 x 24ml
HRP Diluent	Red	1 x 12ml	1 x 23ml
ICAM3 Control	Silver	2 vials	4 vials
ICAM3 Microplate (12 x 8 well strips)		1 unit	2 units
ICAM3 Standard	Yellow	2 vials	4 vials
Stop Reagent	Black	1 x 11ml	2 x 11ml
Streptavidin-HRP		2 x 5µl	4 x 5µl

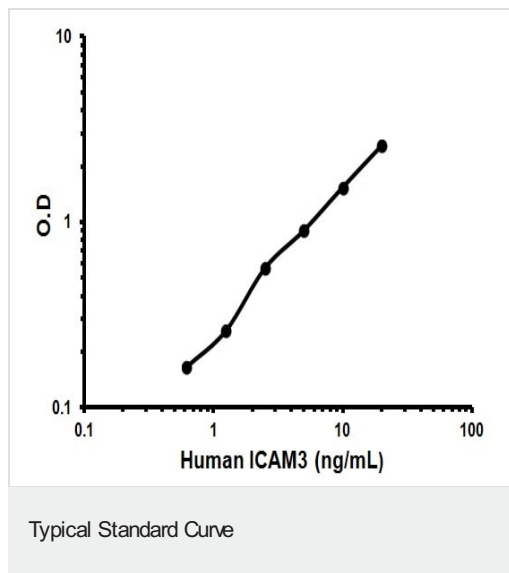
Relevance

ICAM3 is a member of the intercellular adhesion molecule family. All ICAM proteins are type I transmembrane glycoproteins, contain 2-9 immunoglobulin-like C2-type domains, and bind to the leukocyte adhesion LFA-1 protein. ICAM3 is constitutively and abundantly expressed by all leucocytes and may be the most important ligand for LFA-1 in the initiation of the immune response. It functions not only as an adhesion molecule, but also as a potent signalling molecule.

Cellular localization

Cell Membrane; single-pass type I membrane protein

Images



Representative Standard Curve using ab46024

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