

Human ICAM2 ELISA Kit (CD102) ab46080

画像数 1

製品の概要

製品名	Human ICAM2 ELISA Kit (CD102)				
検出方法	Colorimetric				
再現性	Intra-Assay (同時再現性)				
	サンプル	N	平均値	SD	CV%
	A	8	23.62U/ml	0.277	1.17%
	B	8	2.96U/ml	0.86	2.9%
	Inter-Assay (日差再現性)				
	サンプル	N	平均値	SD	CV%
	A	33	23.26U/ml	0.668	2.87%
	B	33	3.12	0.218	6.98%
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	= 0.2 U/ml				
検出範囲	0.75 U/ml - 24 U/ml				
全工程の試験時間	1h 45m				
ステップ	Multiple steps standard assay				
種交差性	交差種: Human				
製品の概要	<p>Abcam's Human ICAM2 <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of soluble Inter Cellular Adhesion Molecule -2 (ICAM2) in Human serum, plasma, buffered solutions or cell culture medium.</p> <p>A monoclonal antibody specific for ICAM2 has been coated onto the wells of the microtiter strips provided. Samples, including standards of known ICAM2 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 ICAM2 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</p>				

colored reaction product. The intensity of this colored product is directly proportional to the concentration of ICAM2 present in the samples.

This kit will recognize both endogenous and recombinant Human ICAM2.

試験プラットフォーム

Microplate

製品の特性

保存方法

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

内容	ラベル	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-ICAM2	Red	1 x 400µl	2 x 400µl
Chromogen TMB Substrate Solution		1 x 11ml	1 x 24ml
Control	Silver	2 vials	4 vials
HRP Diluent	Red	1 x 12ml	1 x 23ml
ICAM2 Microplate (12 x 8 well strips)		1 unit	2 units
ICAM2 Standard (Lyophilized)	Yellow	2 vials	4 vials
Stop Reagent	Black	1 x 11ml	2 x 11ml
Streptavidin-HRP		2 x 5µl	4 x 5µl

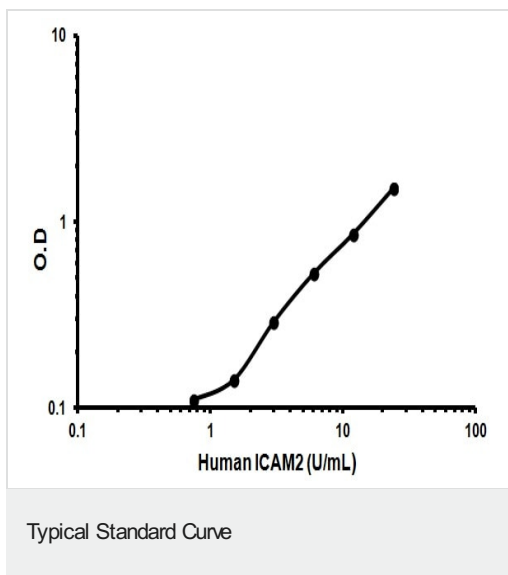
関連性

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.

細胞内局在

Type 1 membrane protein

画像



Representative Standard Curve using ab46080

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