



Product datasheet

Human ICAM3 ELISA Kit (CD50) ab46024

画像数 1

製品の概要

製品名	Human ICAM3 ELISA Kit (CD50)				
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%
	Sample A	8	19.75ng/ml	0.172	0.86%
	Sample B	8	2.83ng/ml	0.099	3.49%
	Inter-Assay (日差再現性)				
	サンプル	N	平均値	SD	CV%
	Sample A	36	19.51ng/ml	0.391	1.99%
	Sample B	36	2.71ng/ml	0.176	6.27%
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	< 0.3 ng/ml				
検出範囲	0.62 ng/ml - 20 ng/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	Abcam's ICAM3 (CD50) Human <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is				

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.

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-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

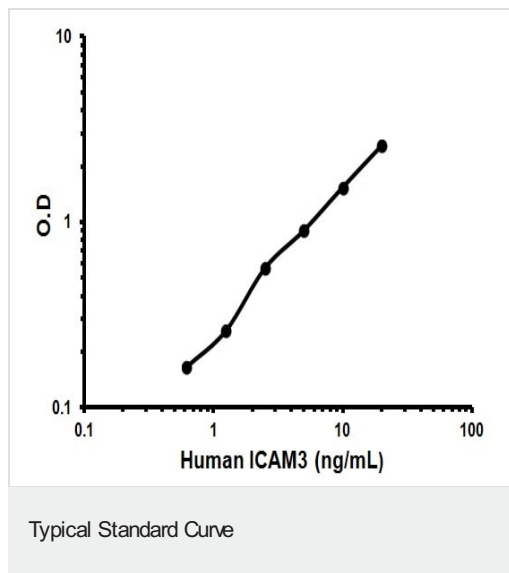
関連性

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.

細胞内局在

Cell Membrane; single-pass type I membrane protein

画像



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.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