

Human ICAM1 ELISA Set (without plates) ab47349

3 References 画像数 1

製品の概要

製品名	Human ICAM1 ELISA Set (without plates)
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	< 0.1 ng/ml
検出範囲	0.25 ng/ml - 8 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	<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>
アプリケーション	適用あり: Sandwich ELISA
試験プラットフォーム	Reagents

製品の特性

保存方法 Store at +4°C. Please refer to protocols.

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

内容	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

機能 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.

配列類似性 Belongs to the immunoglobulin superfamily. ICAM family.
Contains 5 Ig-like C2-type (immunoglobulin-like) domains.

翻訳後修飾 Monoubiquitinated, which is promoted by MARCH9 and leads to endocytosis.

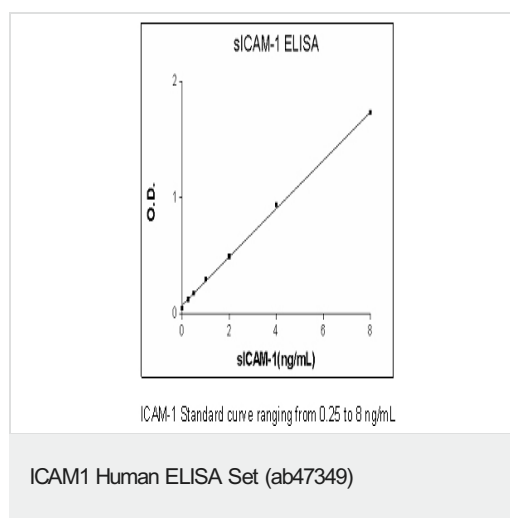
細胞内局在 Membrane.

アプリケーション

The Abpromise guarantee Abpromise保証は、次のテスト済みアプリケーションにおけるab47349の使用に適用されます
アプリケーションノートには、推奨の開始希釈率がありますが、適切な希釈率につきましてはご確認ください。

アプリケーション	Abreviews	特記事項
Sandwich ELISA		Use at an assay dependent concentration.

画像



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