

Human CD86 ELISA Kit ab45921

2 References 画像数 4

製品の概要

製品名	Human CD86 ELISA Kit				
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	37.25U/ml	0.329	0.88%
	Sample B	8	5.09U/ml	0.193	3.79%
	Inter-Assay (日差再現性)				
	サンプル	N	平均値	SD	CV%
	Sample A	39	37.63U/ml	0.611	1.62%
	Sample B	39	4.92U/ml	0.348	7.06%
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	= 0.6 U/ml				
検出範囲	1.187 U/ml - 38 U/ml				
全工程の試験時間	2h 45m				
ステップ	Multiple steps standard assay				
種交差性	交差種: Human				
製品の概要	Abcam's Human CD86 <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for				

colored reaction product. The intensity of this colored product is directly proportional to the concentration of CD86 present in the samples.

This kit will recognize both endogenous and recombinant Human CD86.

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	2 x 96 tests	1 x 96 tests
10X Standard Diluent Buffer	Black	1 x 25ml	1 x 15ml
200X Wash Buffer	White	2 x 10ml	1 x 10ml
Biotinylated Antibody Diluent	Red	1 x 13ml	1 x 7ml
Biotinylated anti-CD86	Red	2 x 0.4ml	1 x 0.4ml
CD86 Microplate (12 x 8 well strips)		2 units	1 unit
CD86 Standard (Lyophilized)	Yellow	4 vials	2 vials
Chromogen TMB Substrate Solution		1 x 24ml	1 x 11ml
HRP Diluent	Red	1 x 23ml	1 x 12ml
Stop Reagent	Black	2 x 11ml	1 x 11ml
Streptavidin-HRP		4 x 5µl	2 x 5µl

機能

Receptor involved in the costimulatory signal essential for T-lymphocyte proliferation and interleukin-2 production, by binding CD28 or CTLA-4. May play a critical role in the early events of T-cell activation and costimulation of naive T-cells, such as deciding between immunity and anergy that is made by T-cells within 24 hours after activation. Isoform 2 interferes with the formation of CD86 clusters, and thus acts as a negative regulator of T-cell activation.

組織特異性

Expressed by activated B-lymphocytes and monocytes.

配列類似性

Contains 1 Ig-like C2-type (immunoglobulin-like) domain.

Contains 1 Ig-like V-type (immunoglobulin-like) domain.

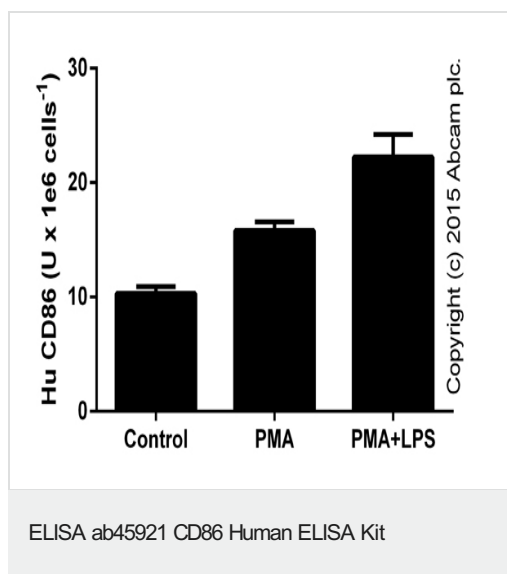
翻訳後修飾

Polyubiquitinated; which is promoted by MARCH8 and results in endocytosis and lysosomal degradation.

細胞内局在

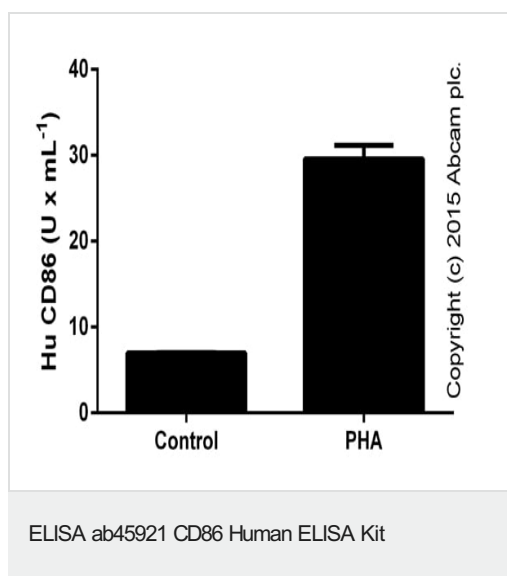
Membrane.

画像

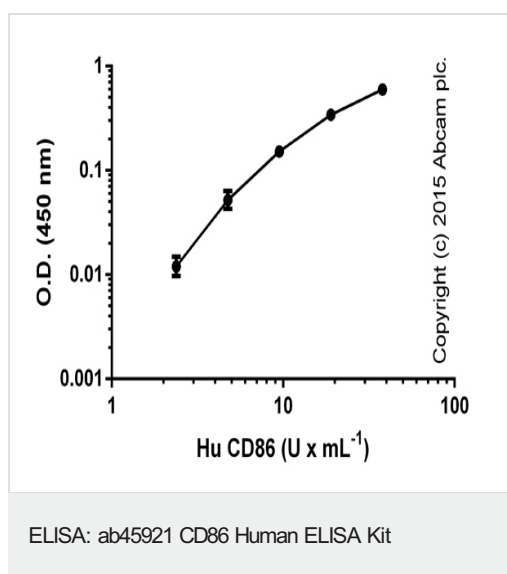


CD86 in U937 lysates from control cells or cells stimulated for 24 hours with 50 ng x mL⁻¹ PMA ([ab120297](#)) or PMA with 1 ug x mL⁻¹ LPS for the last 6 hours (duplicates; +/- SD).

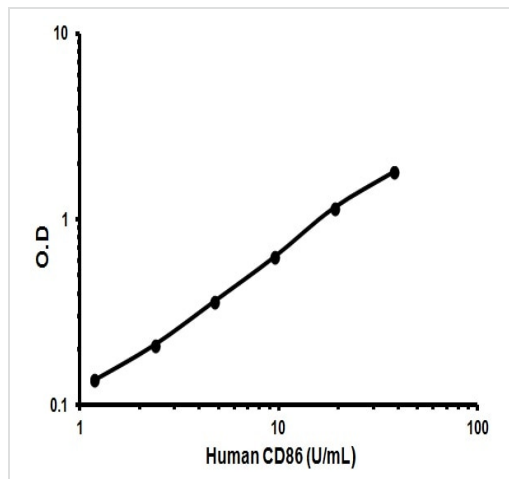
2e6 snap frozen cells were lysed in 0.3 mL RIPA lysis buffer on ice, sonicated in an ice cold water bath for 12 intervals for a total of 3.5 minutes. The supernatants were cleared by centrifugation (max speed, 10 minutes, 4 degrees), and tested undiluted or diluted 1/4.



CD86 measured in undiluted 72 hours culture supernatants from control of treated (5 ug x mL⁻¹ PHA; Sigma) PBMCs (1e6 cells x mL⁻¹; duplicates +/- SD).



Standard curve with background signal subtracted (duplicates; +/- SD).



Representative Standard Curve using ab45921

Typical Standard Curve

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