

Human CD86 ELISA Kit ab45921

2 References 4 图像

概述

产品名称

检测方法

精确度

人CD86 ELISA试剂盒

Colorimetric

批次内

样品	n	Mean	SD	CV%
Sample A	8	37.25U/ml	0.329	0.88%
Sample B	8	5.09U/ml	0.193	3.79%

批次间

样品	n	Mean	SD	CV%
Sample A	39	37.63U/ml	0.611	1.62%
Sample B	39	4.92U/ml	0.348	7.06%

样品类型

检测类型

灵敏度

范围

检测时间

实验步骤

种属反应性

Cell culture supernatant, Serum, Plasma

Sandwich (quantitative)

= 0.6 U/ml

1.187 U/ml - 38 U/ml

2h 45m

Multiple steps standard assay

与反应: Human

产品概述

Abcam's Human CD86 *in vitro* 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.

平台 Microplate

性能

存放说明 Store at +4°C. Please refer to protocols.

组 件	标识符	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

功能 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.

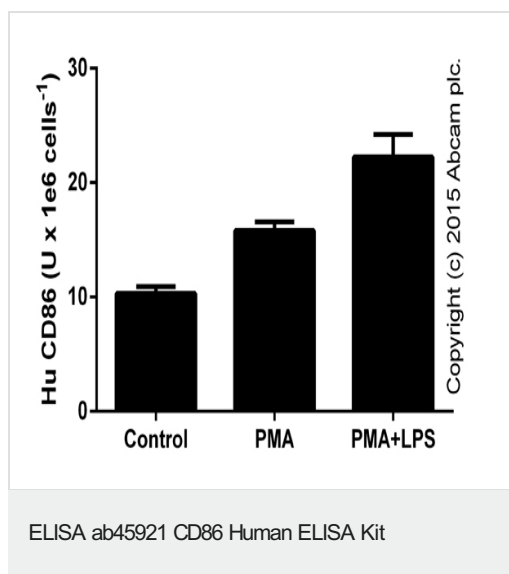
组织特异性 Expressed by activated B-lymphocytes and monocytes.

序列相似性 Contains 1 Ig-like C2-type (immunoglobulin-like) domain.
Contains 1 Ig-like V-type (immunoglobulin-like) domain.

翻译后修饰 Polyubiquitinated; which is promoted by MARCH8 and results in endocytosis and lysosomal degradation.

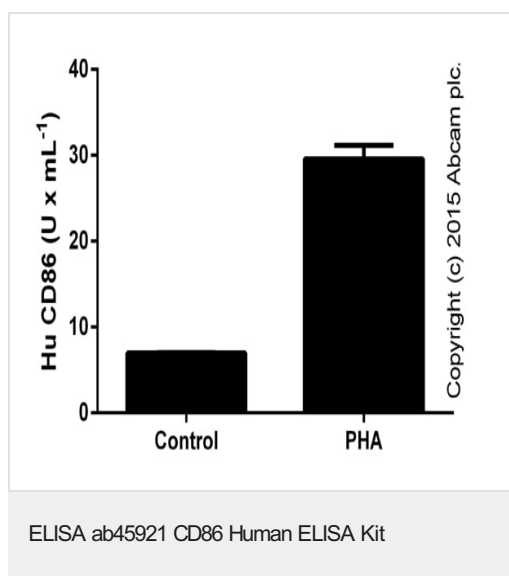
细胞定位 Membrane.

图片

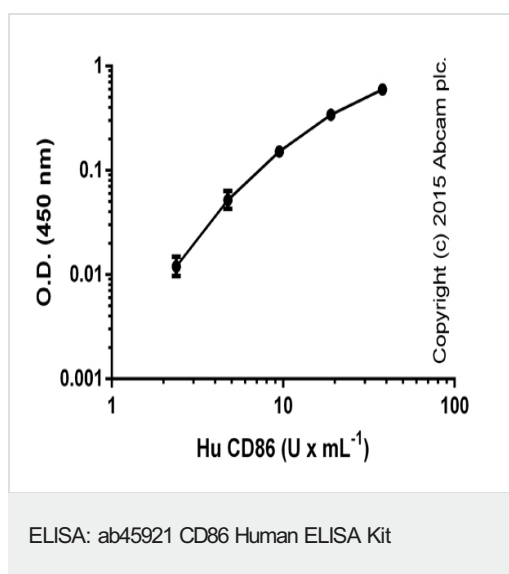


CD86 in U937 lysates from control cells or cells stimulated for 24 hours with 50 ng x mL⁻¹ PMA ([ab120297](#)) or PMA with 1 ug x mL⁻¹ LPS for the last 6 hours (duplicates; +/- SD).

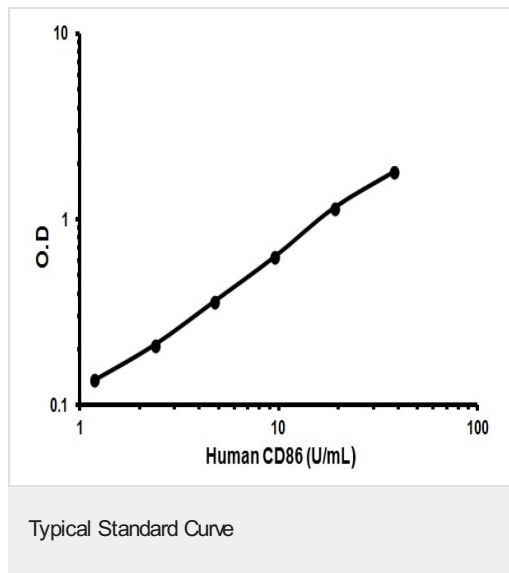
2e6 snap frozen cells were lysed in 0.3 mL RIPA lysis buffer on ice, sonicated in an ice cold water bath for 12 intervals for a total of 3.5 minutes. The supernatants were cleared by centrifugation (max speed, 10 minutes, 4 degrees), and tested undiluted or diluted 1/4.



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

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.cn/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors