

Human IFN gamma High Sensitivity ELISA Kit ab46048

3 References 画像数 5

製品の概要

製品名	Human IFN gamma High Sensitivity 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%
	Overall				3.9%
	Inter-Assay (日差再現性)				
	サンプル	N	平均値	SD	CV%
	Overall				8.6%
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	< 0.69 pg/ml				
検出範囲	0.78 pg/ml - 25 pg/ml				
添加回収試験	特定サンプルでの回収試験				
	サンプルの種類		平均 %	測定範囲	
	Cell culture supernatant		108	% - %	
	Serum		92	% - %	
	Plasma		72	% - %	
全工程の試験時間	3h 00m				
ステップ	Multiple steps standard assay				
種交差性	交差種: Human				
製品の概要	Abcam's IFN gamma Human High Sensitivity <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of IFN gamma in Human sera, plasmas, buffered solutions and cell culture media.				

A monoclonal antibody specific for IFN gamma has been coated onto the wells of the microtiter strips provided. Samples, including standards of known IFN gamma 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 IFN gamma 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 colored reaction product. The intensity of this colored product is directly proportional to the concentration of IFN gamma present in the samples.

This kit will recognize both endogenous and recombinant Human IFN gamma.

Get results in 90 minutes with Human IFN gamma ELISA Kit ([ab174443](#)) from our SimpleStep ELISA® range.

## 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.

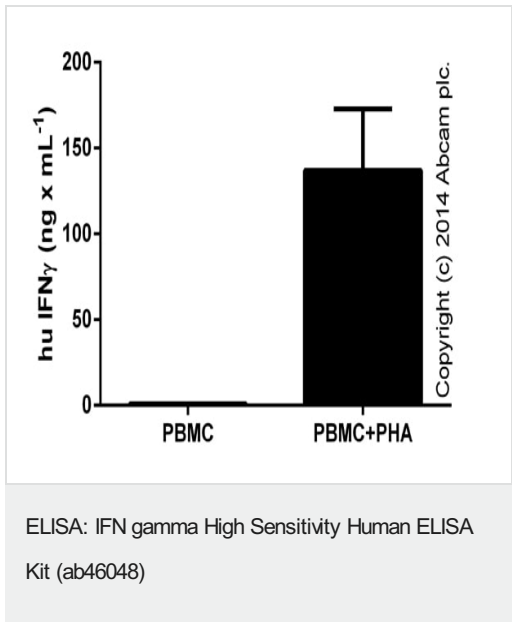
内容	ラベル	2 x 96 tests	1 x 96 tests
10X Standard Diluent Buffer	Black	1 x 25ml	1 x 15ml
200X Wash Buffer	White	2 x 10ml	1 x 10ml
Amplification Diluent	Brown & blue spot	1 x 25ml	1 x 15ml
Amplifier	Yellow	2 x 200µl	1 x 200µl
Biotinylated Antibody Diluent	Red	1 x 13ml	1 x 7.5ml
Biotinylated anti-IFN gamma	Red	2 x 0.4ml	1 x 0.4ml
Chromogen TMB Substrate Solution		1 x 24ml	1 x 11ml
HRP Diluent	Red	2 x 23ml	1 x 23ml
IFN gamma Microplate (12 x 8 well strips)		2 units	1 unit
IFN gamma Standard (Lyophilized)	Yellow	4 vials	2 vials
Plastic Plate Covers		4 units	2 units
Stop Reagent	Black	2 x 11ml	1 x 11ml
Streptavidin-HRP		4 x 5µl	2 x 5µl

## 機能

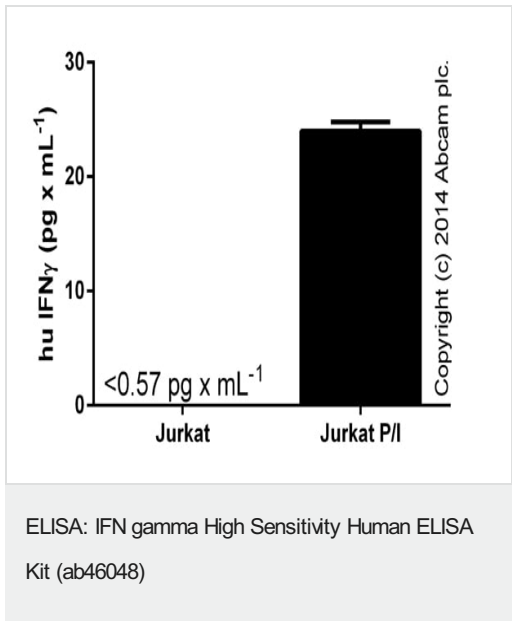
Produced by lymphocytes activated by specific antigens or mitogens. IFN-gamma, in addition to having antiviral activity, has important immunoregulatory functions. It is a potent activator of macrophages, it has antiproliferative effects on transformed cells and it can potentiate the antiviral

	and antitumor effects of the type I interferons.
組織特異性	Released primarily from activated T lymphocytes.
関連疾患	In Caucasians, genetic variation in IFNG is associated with the risk of aplastic anemia (AA) [MIM:609135]. AA is a rare disease in which the reduction of the circulating blood cells results from damage to the stem cell pool in bone marrow. In most patients, the stem cell lesion is caused by an autoimmune attack. T-lymphocytes, activated by an endogenous or exogenous, and most often unknown antigenic stimulus, secrete cytokines, including IFN-gamma, which would in turn be able to suppress hematopoiesis.
配列類似性	Belongs to the type II (or gamma) interferon family.
翻訳後修飾	Proteolytic processing produces C-terminal heterogeneity, with proteins ending alternatively at Gly-150, Met-157 or Gly-161.
細胞内局在	Secreted.

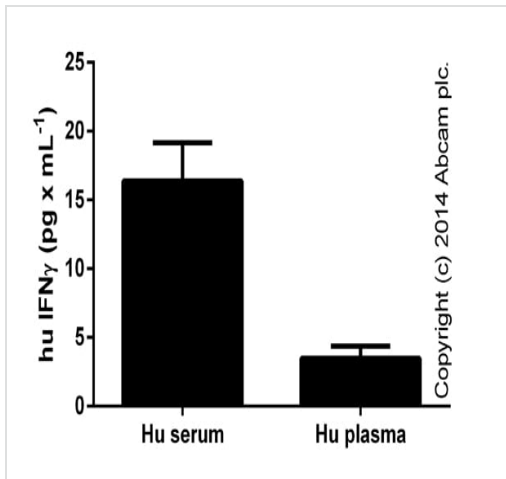
画像



IFN gamma measured in cell culture supernatants from control or treated (48 hours; 2% PHA-M; LifeTechnologies) PBMCs, diluted 1/100-1/10000 (duplicates +/- SD).

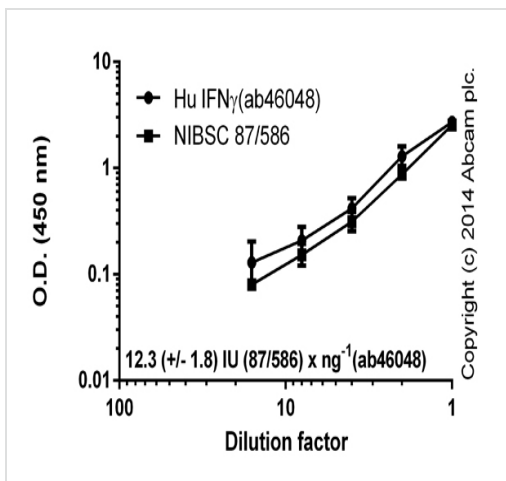


IFN gamma measured in undiluted cell culture supernatants from control or stimulated (48 hours with 50 ng x mL<sup>-1</sup> PMA ([ab120297](#)) and 1  $\mu$ M ionomycin ([ab120116](#))) Jurkat cells. Control sample was below limit of detection (<0.57 pg x mL<sup>-1</sup>) (duplicates +/- SD).



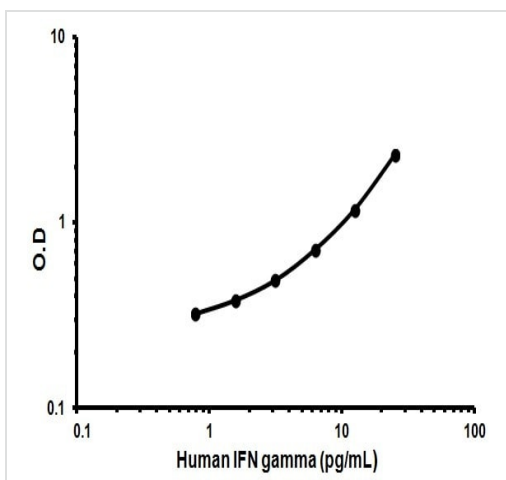
ELISA: IFN gamma High Sensitivity Human ELISA  
Kit (ab46048)

IFN gamma measured in undiluted human biological fluids; mouse and rat samples (plasma and serum) were below limit of detection (duplicates  $\pm$  SD).



ELISA: IFN gamma High Sensitivity Human ELISA  
Kit (ab46048)

Dilution curves of IFN gamma (ab46048) and NIBSC standard (87/586). One ng of standard IFN gamma corresponds to 12.3 ( $\pm$  1.3) IU NIBSC 87/586. Background signal subtracted (duplicates;  $\pm$  SD).



Typical Standard Curve

Representative Standard Curve using ab46048

**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