

Human Interferon gamma ELISA Kit (IFNG) ab100537

3 References 画像数 2

製品の概要

製品名	Human Interferon gamma ELISA Kit (IFNG)
検出方法	Colorimetric
サンプルの種類	Cell culture supernatant, Plasma
アッセイタイプ	Sandwich (quantitative)
検出感度	< 15 pg/ml
検出範囲	20.6 pg/ml - 15000 pg/ml
添加回収試験	> 87 %

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	94.53	84% - 104%
Plasma	86.82	81% - 102%

ステップ	Multiple steps standard assay
種交差性	交差種: Human
製品の概要	

Abcam’s Interferon gamma (IFNG) Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Human Interferon gamma- in plasma, and cell culture supernatants. (Human Interferon gamma concentration is pretty low in normal plasma, it may not be detected in this assay).

This assay employs an antibody specific for Human Interferon gamma coated on a 96-well plate. Standards and samples are pipetted into the wells and Interferon gamma present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human Interferon gamma antibody is added. After washing away unbound biotinylated antibody, HRP-conjugated streptavidin is pipetted to the wells. The wells are again washed, a TMB substrate solution is added to the wells and color develops in proportion to the amount of Interferon gamma bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

We have not been able to detect endogenous Human Interferon gamma in normal serum with ab100537, only in serum spiked with Human Interferon gamma.

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

特記事項

Optimization may be required with urine samples

試験プラットフォーム

Microplate

製品の特性

保存方法

Store at -20°C. Please refer to protocols.

内容	1 x 96 tests
20X Wash Buffer	1 x 25ml
400X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent B	1 x 15ml
Assay Diluent A	1 x 30ml
Biotinylated anti-Human Interferon gamma	2 vials
Interferon gamma Microplate (12 x 8 wells)	1 unit
Recombinant Human Interferon gamma Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

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.

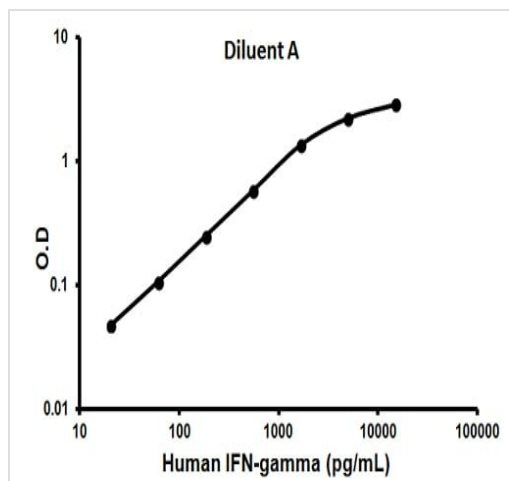
翻訳後修飾

Proteolytic processing produces C-terminal heterogeneity, with proteins ending alternatively at Gly-150, Met-157 or Gly-161.

細胞内局在

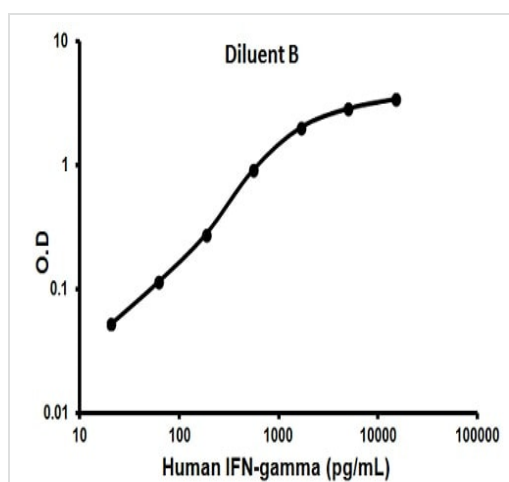
Secreted.

画像



Representative standard curve using ab100537

Typical standard curve



Representative standard curve using ab100537

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