

Rat IFN gamma ELISA Kit ab113349

3 References [画像数 2](#)

製品の概要

製品名	Rat IFN gamma ELISA Kit
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	< 100 pg/ml
検出範囲	123 pg/ml - 30000 pg/ml
添加回収試験	98 %

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	99.29	89% - 111%
Plasma	92.63	78% - 104%

ステップ Multiple steps standard assay

種交差性 交差種: Rat

製品の概要

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

This assay employs an antibody specific for IFN gamma coated on a 96-well plate. Standards and samples are pipetted into the wells and IFN gamma present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Rat IFN 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 IFN 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 the endogenous Rat IFN-gamma in normal serum with ab113349, only in serum spiked with Rat IFN-gamma.

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
200X HRP-Streptavidin Concentrate	1 x 200µl
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
Assay Diluent A	1 x 30ml
Biotinylated anti-Rat IFN Gamma	2 vials
IFN gamma Microplate (12 x 8 wells)	1 x 96 units
Recombinant rat IFN 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.

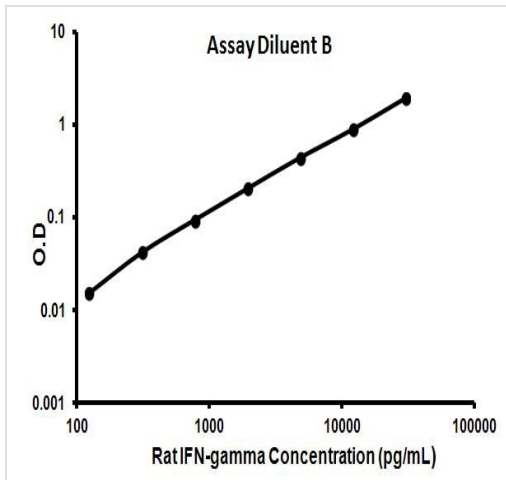
翻訳後修飾

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

細胞内局在

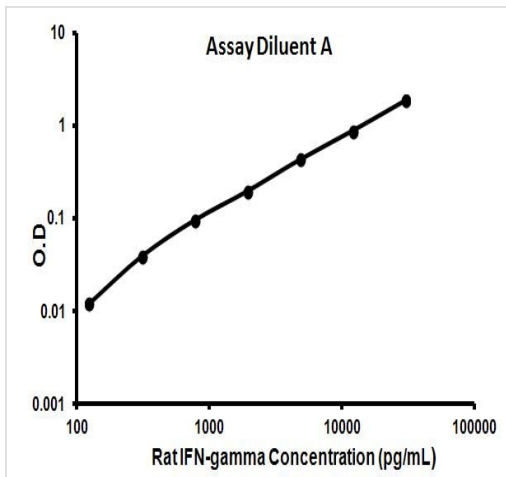
Secreted.

画像



Representative standard curve using ab113349

Typical data B



Representative standard curve using ab113349

Typical data A

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