

# Rat IFN gamma ELISA Kit ab113349

3 References 2 图像

概述

产品名称	大鼠IFN gamma ELISA试剂盒
检测方法	Colorimetric
样品类型	Cell culture supernatant, Plasma
检测类型	Sandwich (quantitative)
灵敏度	< 100 pg/ml
范围	123 pg/ml - 30000 pg/ml
回收率	98 %

特定样本回收率

样品类型	平均%	范围
Cell culture supernatant	99.29	89% - 111%
Plasma	92.63	78% - 104%

实验步骤 Multiple steps standard assay

种属反应性 与反应: Rat

产品概述 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.

平台 Microplate

性能

存放说明 Store at -20°C. Please refer to protocols.

组件	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

功能 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.

组织特异性 Released primarily from activated T lymphocytes.

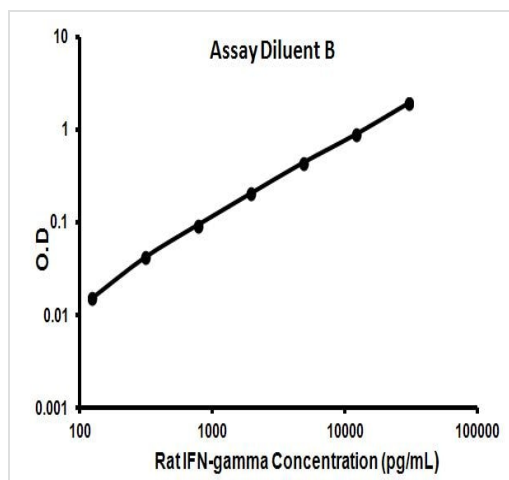
疾病相关 In Caucasians, genetic variation in IFNG is associated with the risk of aplastic anemia (AA) [MIM:609135]. AA is a rare disease in which the reduction of the circulating blood cells results from damage to the stem cell pool in bone marrow. In most patients, the stem cell lesion is caused by an autoimmune 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.

序列相似性 Belongs to the type II (or gamma) interferon family.

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

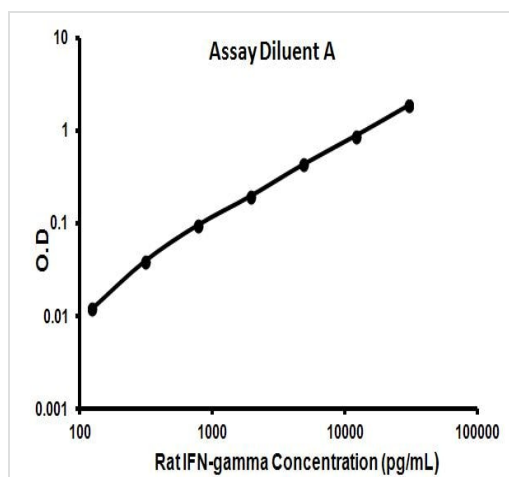
细胞定位 Secreted.

图片



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.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