

### Mouse Interferon gamma ELISA Kit ab46081

★★★★★ [1 Abreviews](#) [5 References](#) [1 图像](#)

#### 概述

**产品名称** 小鼠Interferon gamma ELISA试剂盒

**检测方法** Colorimetric

**精确度** 批次内

样品	n	Mean	SD	CV%
A	20	315.4pg/ml	14.5	4.6%
B	19	588.6pg/ml	19.1	4.6%

批次间

样品	n	Mean	SD	CV%
A	9	128.8pg/ml	7.4	5.7%
B	10	546.9pg/ml	25.2	4.6%

**样品类型** Milk, Serum

**检测类型** Sandwich (quantitative)

**灵敏度** < 15 pg/ml

**范围** 31.25 pg/ml - 1000 pg/ml

**检测时间** 3h 45m

**实验步骤** Multiple steps standard assay

**种属反应性** **与反应:** Mouse

**产品概述** Mouse interferon gamma ELISA kit (IFN Gamma) is designed for the quantitative measurement of IFN gamma in mouse serum, buffered solutions and cell culture medium.

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 mouse IFN gamma.

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

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 7.5ml
Chromogen TMB Substrate Solution		1 x 24ml	1 x 11ml
HRP Diluent	Red	1 x 23ml	1 x 12ml
IFN gamma Microplate (12 x 8 well strips)		2 units	1 unit
IFN gamma Standard (Lyophilized)	Yellow	4 vials	2 vials
IFN $\gamma$ Biotinylated detection antibody		2 x 400 $\mu$ l	1 x 400 $\mu$ l
Stop Reagent	Black	2 x 11ml	1 x 11ml
Streptavidin-HRP		4 x 5 $\mu$ l	2 x 5 $\mu$ l

**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.

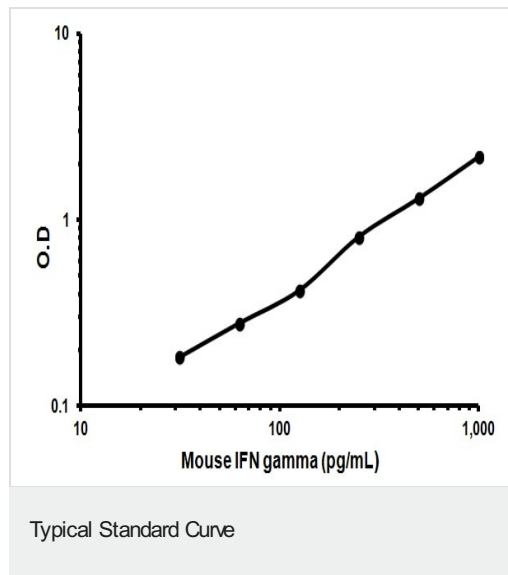
**组织特异性** Released primarily from activated T lymphocytes.

**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.

## 图片



Representative Standard Curve using ab46081

**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