

### Mouse Interferon gamma ELISA Kit (IFNG) ab100689

★★★★☆ [2 Abreviews](#) [28 References](#) [2 Images](#)

#### Overview

<b>Product name</b>	Mouse Interferon gamma ELISA Kit (IFNG)
<b>Detection method</b>	Colorimetric
<b>Sample type</b>	Cell culture supernatant, Serum, Plasma
<b>Assay type</b>	Sandwich (quantitative)
<b>Sensitivity</b>	= 1.3 pg/ml
<b>Range</b>	2.74 pg/ml - 2000 pg/ml
<b>Recovery</b>	94 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	94.75	83% - 103%
Serum	95.38	83% - 103%
Plasma	93.49	82% - 102%

**Assay duration** Multiple steps standard assay

**Species reactivity** **Reacts with:** Mouse

**Product overview** Abcam's Interferon Gamma Mouse ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Mouse Interferon Gamma in serum, plasma and cell culture supernatants.

This assay employs an antibody specific for Mouse 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-Mouse 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.

**Platform** Microplate

## Properties

### Storage instructions

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

Components	1 x 96 tests
20X Wash Buffer	1 x 25ml
440X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent B	1 x 15ml
Assay Diluent A	1 x 30ml
Biotinylated anti-Mouse Interferon gamma	2 vials
Interferon gamma Microplate (12 strips x 8 wells)	1 unit
Recombinant Mouse Interferon gamma Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

### Function

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.

### Tissue specificity

Released primarily from activated T lymphocytes.

### Involvement in disease

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.

### Sequence similarities

Belongs to the type II (or gamma) interferon family.

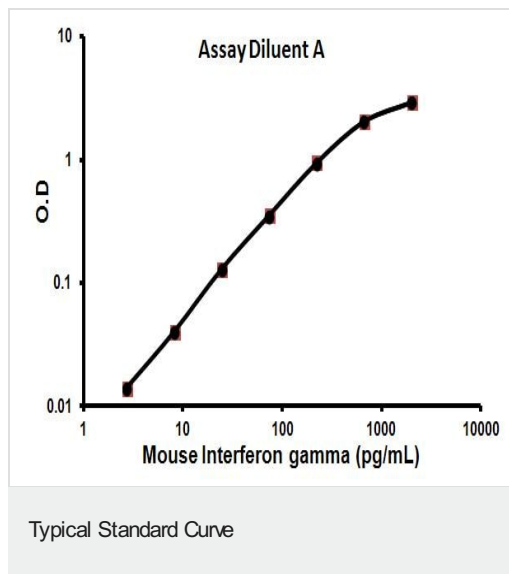
### Post-translational modifications

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

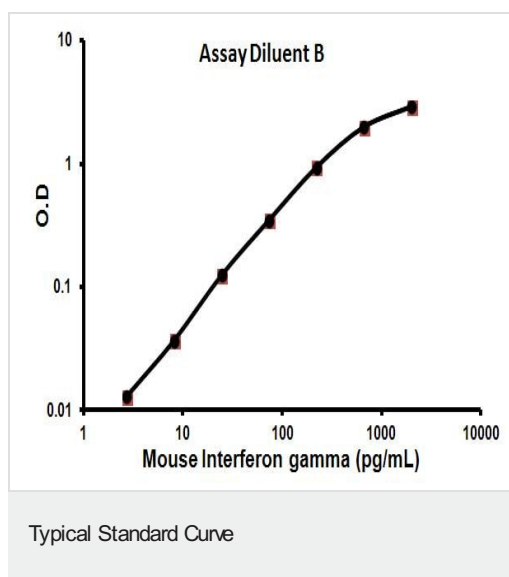
### Cellular localization

Secreted.

## Images



Representative Standard Curve using ab100689



Representative Standard Curve using ab100689

**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.com/abpromise> or contact our technical team.

#### **Terms and conditions**

---

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors