

Human Interferon gamma ELISA Kit (IFNG) ab108863

1 Image

Overview

Product name	Human Interferon gamma ELISA Kit (IFNG)			
Detection method	Colorimetric			
Precision	Intra-assay			
	Sample	n	Mean	SD
	Overall			5.2%
	Inter-assay			
	Sample	n	Mean	SD
	Overall			9.8%
Sample type	Cell culture supernatant, Serum, Plasma			
Assay type	Sandwich (quantitative)			
Sensitivity	> 0.01 ng/ml			
Range	0.016 ng/ml - 1 ng/ml			
Recovery	101 %			
Assay time	5h 00m			
Assay duration	Multiple steps standard assay			
Species reactivity	Reacts with: Human			
Product overview	Abcam's Interferon gamma (IFNG) Human <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of Interferon gamma levels in cell culture supernatants, serum and plasma.			

An Interferon gamma specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently an Interferon gamma specific biotinylated detection antibody is added and then followed by washing with wash buffer. Streptavidin-Peroxidase Conjugate is added and unbound conjugates are washed away with wash buffer. TMB is then used to visualize Streptavidin-Peroxidase enzymatic reaction. TMB is catalyzed by Streptavidin-Peroxidase to produce a blue color product that changes into yellow after adding acidic stop solution. The density of yellow coloration is directly proportional to the amount of Interferon gamma captured in plate.

The entire kit may be stored at -20°C for long term storage before reconstitution - Avoid repeated freeze-thaw cycles.

Platform Microplate

Properties

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

Components	1 x 96 tests
100X Streptavidin-Peroxidase Conjugate	1 x 80µl
10X Diluent M Concentrate	1 x 20ml
1X Standard Diluent	1 x 2ml
20X Wash Buffer Concentrate	2 x 30ml
40X Biotinylated Human Interferon gamma Antibody	1 x 150µl
Chromogen Substrate	1 x 8ml
Interferon gamma Microplate (12 x 8 well strips)	1 unit
Interferon gamma Standard	1 vial
Sealing Tapes	3 units
Stop Solution	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.

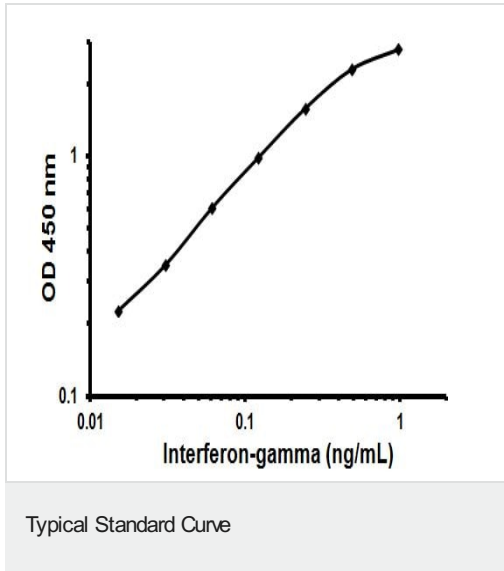
Involvement in disease In Caucasians, genetic variation in IFNG is associated with the risk of aplastic anemia (AA) [MIM:609135]. AA is a rare disease in which the reduction of the circulating blood cells results from damage to the stem cell pool in bone marrow. In most patients, the stem cell lesion 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 ab108863.

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