

Mouse Interferon gamma ELISA Set (without plates) ab47618

[1 Image](#)

Overview

Product name	Mouse Interferon gamma ELISA Set (without plates)
Sample type	Cell culture supernatant, Serum, Plasma
Assay type	Sandwich (quantitative)
Assay duration	Multiple steps standard assay
Species reactivity	Reacts with: Mouse
Product overview	<p>Abcam's IFNγ Murine <i>in vitro</i> ELISA Set is designed for the quantitative measurement of Murine IFNγ in cell culture supernatants, buffered solutions, serum, plasma and other body fluids.</p> <p>The murine IFN-gamma ELISA Set includes both the capture and the detection antibodies as well as Horse-Radish peroxidase conjugated Streptavidin (HRP-Strep) and TMB which are the reagents used in the colour development step.</p> <p>This ELISA Set is intended to measure murine IFN-gamma in cell -culture supernatants, as well as in serum samples. The enclosed procedures provide general guidelines for ELISA preparation. This procedure may need some optimization depending on the plastic, matrix, standard range.</p> <p>Recognizes both natural and recombinant Murine Interferon gamma Range: 31.25 pg/ml - 1000 pg/ml Sensitivity < 15 pg/ml.</p>
Tested applications	Suitable for: Sandwich ELISA
Platform	Reagents

Properties

Storage instructions Store at +4°C. Please refer to protocols.

Components	Identifier	15 x 96 tests	5 x 96 tests	10 x 96 tests
Chromogen TMB Substrate Solution		6 vials	2 vials	4 vials
Detection biotinylated anti-Murine IFN gamma antibody		3 vials	1 vial	2 vials

Components	Identifier	15 x 96 tests	5 x 96 tests	10 x 96 tests
IFN gamma Standard (Lyophilized)	Yellow	15 vials	5 vials	10 vials
Murine IFN gamma Capture Antibody		3 x 0.25ml	1 x 0.25ml	2 x 0.25ml
Streptavidin-HRP		3 vials	1 vial	2 vials

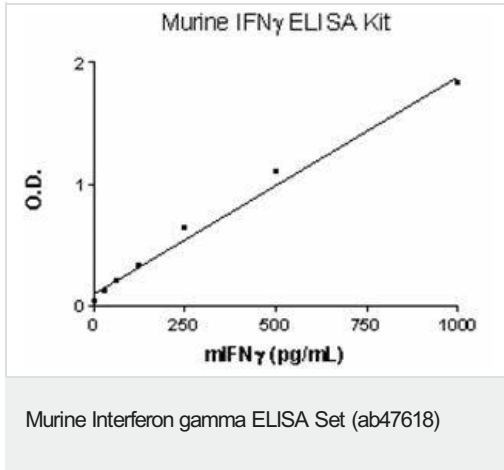
Function	Produced by lymphocytes activated by specific antigens or mitogens. IFN-gamma, in addition to having antiviral activity, has important immunoregulatory functions. It is a potent activator of macrophages, it has antiproliferative effects on transformed cells and it can potentiate the antiviral 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.

Applications

The Abpromise guarantee Our **Abpromise guarantee** covers the use of ab47618 in the following tested applications. The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

Application	Abreviews	Notes
Sandwich ELISA		Use at an assay dependent concentration.

Images



Murine IFN-gamma Standard curve ranging from 31.25 to 1000 pg/mL

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