

Mouse Interferon gamma ELISA Set (without plates) ab47618

画像数 1

製品の概要

製品名	Mouse Interferon gamma ELISA Set (without plates)
サンプルの種類	Cell culture supernatant, Serum, Plasma
アッセイタイプ	Sandwich (quantitative)
ステップ	Multiple steps standard assay
種交差性	交差種: Mouse
製品の概要	<p>Abcam's IFN<math>\gamma</math> Murine <i>in vitro</i> ELISA Set is designed for the quantitative measurement of Murine IFN<math>\gamma</math> 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 <b>Range:</b> 31.25 pg/ml - 1000 pg/ml <b>Sensitivity</b> &lt; 15 pg/ml.</p>
アプリケーション	適用あり: Sandwich ELISA
試験プラットフォーム	Reagents

製品の特性

保存方法 Store at +4°C. Please refer to protocols.

内容	ラベル	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
IFN gamma Standard (Lyophilized)	Yellow	15 vials	5 vials	10 vials

内容	ラベル	15 x 96 tests	5 x 96 tests	10 x 96 tests
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

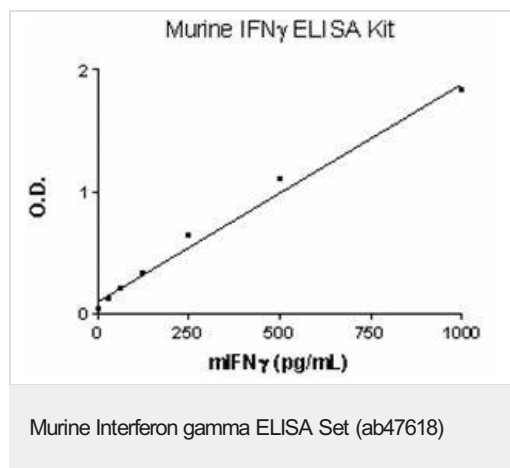
<b>機能</b>	Produced by lymphocytes activated by specific antigens or mitogens. IFN-gamma, in addition to having antiviral activity, has important immunoregulatory functions. It is a potent activator of macrophages, it has antiproliferative effects on transformed cells and it can potentiate the antiviral and antitumor effects of the type I interferons.
<b>組織特異性</b>	Released primarily from activated T lymphocytes.
<b>関連疾患</b>	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.
<b>配列類似性</b>	Belongs to the type II (or gamma) interferon family.
<b>翻訳後修飾</b>	Proteolytic processing produces C-terminal heterogeneity, with proteins ending alternatively at Gly-150, Met-157 or Gly-161.
<b>細胞内局在</b>	Secreted.

## アプリケーション

**The Abpromise guarantee**      **Abpromise保証は、次のテスト済みアプリケーションにおけるab47618の使用に適用されます**  
 アプリケーションノートには、推奨の開始希釈率がありますが、適切な希釈率につきましてはご検討ください。

アプリケーション	Abreviews	特記事項
<b>Sandwich ELISA</b>		Use at an assay dependent concentration.

## 画像



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.co.jp/abpromise> or contact our technical team.

### **Terms and conditions**

---

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