

## Monkey IFN gamma ELISA Kit ab270895

SimpleStep ELISA

画像数 4

### 製品の概要

**製品名** Monkey IFN gamma ELISA Kit

**検出方法** Colorimetric

**再現性** Intra-Assay (同時再現性)

サンプル	N	平均値	SD	CV%
Supernatant	8			5%

Inter-Assay (日差再現性)

サンプル	N	平均値	SD	CV%
Supernatant	3			12%

**サンプルの種類** Cell culture supernatant, Serum, Hep Plasma, EDTA Plasma, Cit plasma, Cerebral Spinal Fluid

**アッセイタイプ** Sandwich (quantitative)

**検出感度** 65 pg/ml

**検出範囲** 0.16 ng/ml - 10 ng/ml

**添加回収試験** 特定サンプルでの回収試験

サンプルの種類	平均 %	測定範囲
Cell culture supernatant	109	97% - 117%
Serum	85	75% - 92%
Hep Plasma	99	80% - 118%
EDTA Plasma	103	95% - 111%
Cit plasma	85	82% - 89%
Cerebral Spinal Fluid	102	98% - 109%

<b>全工程の試験時間</b>	1h 30m
<b>ステップ</b>	One step assay
<b>種交差性</b>	<b>交差種:</b> Monkey
<b>製品の概要</b>	<p>Monkey IFN ELISA kit (ab270895) is a single-wash 90 min sandwich ELISA designed for the quantitative measurement of IFN protein in monkey serum, plasma and cell culture supernatant. It uses our proprietary SimpleStep ELISA® technology. Quantitate monkey IFN with 65 pg/mL sensitivity.</p> <p>SimpleStep ELISA® technology employs capture antibodies conjugated to an affinity tag that is recognized by the monoclonal antibody used to coat our SimpleStep ELISA® plates. This approach to sandwich ELISA allows the formation of the antibody-analyte sandwich complex in a single step, significantly reducing assay time. See the SimpleStep ELISA® protocol summary in the image section for further details. Our SimpleStep ELISA® technology provides several benefits:</p> <ul style="list-style-type: none"> <li>-Single-wash protocol reduces assay time to 90 minutes or less</li> <li>-High sensitivity, specificity and reproducibility from superior antibodies</li> <li>-Fully validated in biological samples</li> <li>-96-wells plate breakable into 12 x 8 wells strips</li> </ul> <p>384-well SimpleStep ELISA® microplate (<b>ab203359</b>) is available to use as an alternative to the 96-well microplate provided with SimpleStep ELISA® kits.</p>
<b>特記事項</b>	<p>IFN gamma (Interferon-gamma), is an important immunoregulatory cytokine in addition of having antiviral activity. 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. IFN gamma is produced by lymphocytes activated by specific antigens or mitogens.</p>
<b>試験プラットフォーム</b>	Pre-coated microplate (12 x 8 well strips)

## 製品の特性

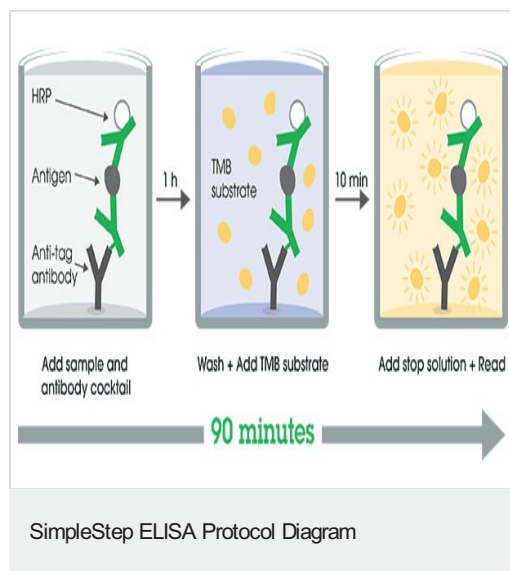
**保存方法** Store at +4°C. Please refer to protocols.

内容	1 x 96 tests
10X Monkey IFN gamma Capture Antibody	1 x 600µl
10X Monkey IFN gamma Detector Antibody	1 x 600µl
10X Wash Buffer PT (ab206977)	1 x 20ml
Antibody Diluent 4BI	1 x 6ml
Monkey IFN gamma Lyophilized Recombinant Protein	2 vials
Plate Seals	1 unit
Sample Diluent 75BP	1 x 20ml
Sample Diluent NS (ab193972)	1 x 12ml

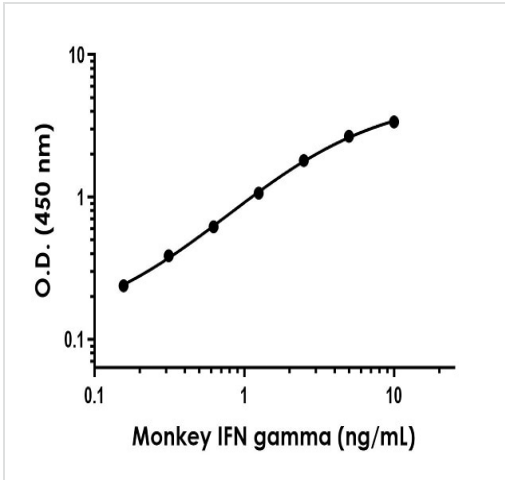
内容	1 x 96 tests
SimpleStep Pre-Coated 96-Well Microplate (ab206978)	1 unit
Stop Solution	1 x 12ml
TMB Development Solution	1 x 12ml

<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.

## 画像

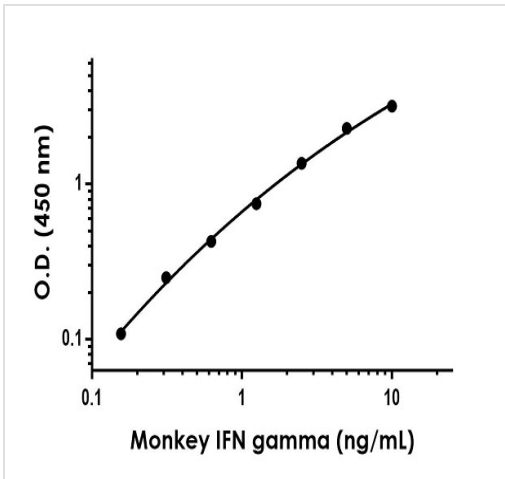


SimpleStep ELISA technology allows the formation of the antibody-antigen complex in one single step, reducing assay time to 90 minutes. Add samples or standards and antibody mix to wells all at once, incubate, wash, and add your final substrate. See protocol for a detailed step-by-step guide.



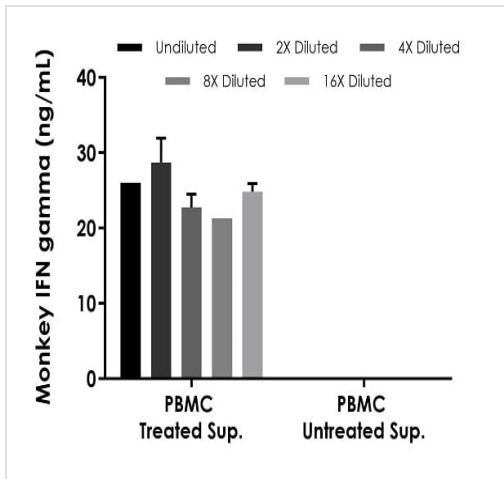
Example of monkey IFN gamma standard curve. Background-subtracted data values (mean +/- SD) are graphed.

Example of monkey IFN gamma standard curve in Sample Diluent NS.



The IFN gamma standard curve was prepared as described in Section 10. Raw data values are shown in the table. Background-subtracted data values (mean +/- SD) are graphed.

Example of monkey IFN gamma standard curve in Sample Diluent 75BP.



Interpolated concentrations of native IFN gamma in cynomolgus monkey PBMC cell culture supernatant samples treated with or without PMA (50 ng/mL) and Ionomycin (1 µg/mL) for 24 hours.

The concentrations of IFN gamma were measured in duplicates, interpolated from the IFN gamma standard curves and corrected for sample dilution. Undiluted samples are as follows: PBMC treated supernatant 10%, and PBMC untreated supernatant 25%. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean IFN gamma concentration was determined to be 25 ng/mL in neat PBMC treated supernatant and undetectable in PBMC untreated supernatant.

**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