

### Human Interferon Alpha 2 ELISA Kit ab233622

リコンビナント SimpleStep ELISA

画像数 5

#### 製品の概要

製品名 Human Interferon Alpha 2 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%
Spiked serum	3			2.7%

Inter-Assay (日差再現性)

サンプル	N	平均値	SD	CV%
Spiked serum	8			9.3%

サンプルの種類 Serum, Hep Plasma, EDTA Plasma, Cit plasma

アッセイタイプ Sandwich (quantitative)

検出感度 56.04 pg/ml

検出範囲 187.5 pg/ml - 12000 pg/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
Serum	103	99% - 109%
Cell culture media	106	100% - 116%
Hep Plasma	104	94% - 108%
EDTA Plasma	87	80% - 93%
Cit plasma	103	102% - 103%

全工程の試験時間 1h 30m

ステップ One step assay

**種交差性**

交差種: Human

**製品の概要**

Human Interferon Alpha 2 ELISA Kit (ab233622) is a single-wash 90 min sandwich ELISA designed for the quantitative measurement of Interferon Alpha 2 protein in cit plasma, edta plasma, hep plasma, and serum. It uses our proprietary SimpleStep ELISA® technology. Quantitate Human Interferon Alpha 2 with 56.04 pg/ml sensitivity.

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:

- Single-wash protocol reduces assay time to 90 minutes or less
- High sensitivity, specificity and reproducibility from superior antibodies
- Fully validated in biological samples
- 96-wells plate breakable into 12 x 8 wells strips

A 384-well SimpleStep ELISA® microplate (**ab203359**) is available to use as an alternative to the 96-well microplate provided with SimpleStep ELISA® kits.

**試験プラットフォーム**

Microplate (12 x 8 well strips)

**製品の特徴****保存方法**

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

内容	1 x 96 tests	10 x 96 tests
10X Human Interferon Alpha 2 Capture Antibody	1 x 600µl	1 x 6000µl
10X Human Interferon Alpha 2 Detector Antibody	1 x 600µl	1 x 6000µl
10X Wash Buffer PT (ab206977)	1 x 20ml	1 x 200ml
Antibody Diluent 5BI	1 x 6ml	10 x 6ml
Human Interferon Alpha 2 Lyophilized Purified Protein	2 vials	2 x 10 vials
Plate Seals	1 unit	1 x 10 units
Sample Diluent NS (ab193972)	1 x 50ml	2 x 250ml
SimpleStep Pre-Coated 96-Well Microplate (ab206978)	1 unit	1 x 10 units
Stop Solution	1 x 12ml	1 x 120ml
TMB Development Solution	1 x 12ml	1 x 120ml

**機能**

Produced by macrophages, IFN-alpha have antiviral activities.

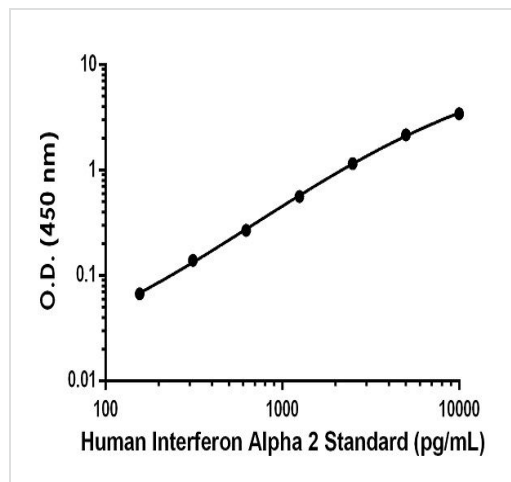
配列類似性

Belongs to the alpha/beta interferon family.

細胞内局在

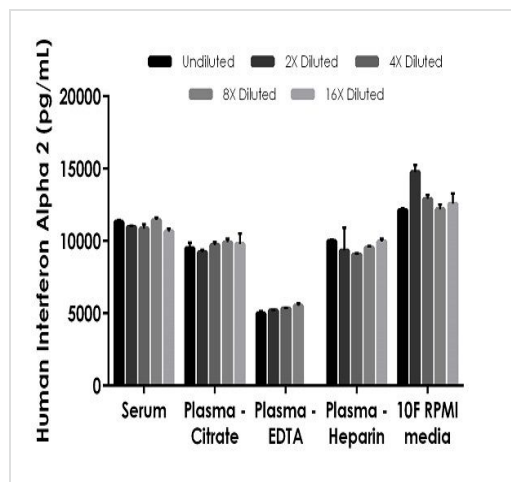
Secreted.

画像



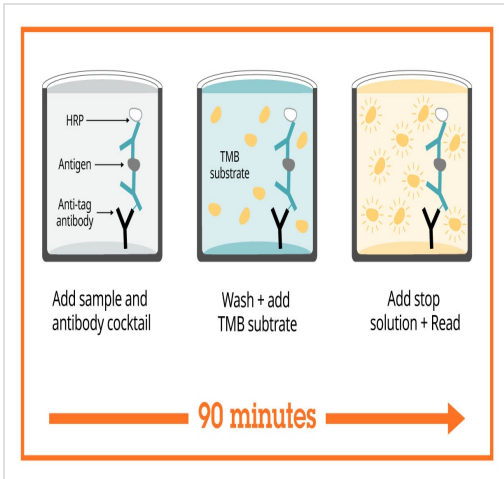
Raw data values are shown in the table. Background-subtracted data values (mean +/- SD) are graphed.

Example of human Interferon Alpha 2 standard curve in Sample Diluent NS



The concentrations of Interferon Alpha 2 were measured in duplicates, interpolated from the Interferon Alpha 2 standard curves.

Interpolated concentrations of spiked Interferon Alpha 2 in human serum, plasma and culture media samples



Sandwich ELISA - Human Interferon Alpha 2 ELISA Kit (ab233622)

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.

Powered by recombinant antibodies

- Research with confidence**  
Consistent and reproducible results
- Long-term and scalable supply**  
Recombinant technology
- Success from the first experiment**  
Confirmed specificity
- Ethical standards compliant**  
Animal-free production

Sandwich ELISA - Human Interferon Alpha 2 ELISA Kit (ab233622)

To learn more about the advantages of recombinant antibodies see [here](#).

**Get more done with SimpleStep ELISA**



**Easy to use**  
Single-wash 90-minute protocol



**Flexible**  
Matched antibody pairs available



**Precision antibodies**  
High sensitivity, specificity and reproducibility



**Scalable**  
Now in 10-pack and 384-well formats

Sandwich ELISA - Human Interferon Alpha 2 ELISA Kit (ab233622)

To learn more about the advantages of SimpleStep ELISA<sup>®</sup> kits see [here](#).

**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