

Mouse IFN beta ELISA Kit ab252363

重组 SimpleStep ELISA

10 References 4 图像

概述

产品名称 小鼠IFN beta ELISA试剂盒

检测方法 Colorimetric

精确度					批次内
	样品	n	Mean	SD	CV%
	Supernatant	8			4.9%

					批次间
	样品	n	Mean	SD	CV%
	Supernatant	3			3.6%

样品类型 Cell culture supernatant, Serum, Cell culture media, EDTA Plasma

检测类型 Sandwich (quantitative)

灵敏度 28 pg/ml

范围 39.1 pg/ml - 2500 pg/ml

回收率			特定样本回收率
	样品类型	平均%	范围
	Cell culture supernatant	97	94% - 100%
	Serum	86	84% - 87%
	Cell culture media	88	78% - 98%
	EDTA Plasma	84	81% - 87%

检测时间 1h 30m

实验步骤 One step assay

种属反应性 与反应: Mouse

产品概述

Mouse IFN beta ELISA Kit (ab252363) is a single-wash 90 min sandwich ELISA designed for the quantitative measurement of IFN beta protein in cell culture supernatant, EDTA plasma, serum, and cell culture media. It uses our proprietary SimpleStep ELISA® technology. Quantitate Mouse IFN beta with 28 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.

平台

Pre-coated microplate (12 x 8 well strips)

性能

存放说明

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

组件	1 x 96 tests
10X Mouse IFN beta (Interferon beta 1) Capture Antibody	1 x 600µl
10X Mouse IFN beta (Interferon beta 1) Detector Antibody	1 x 600µl
10X Wash Buffer PT (ab206977)	1 x 20ml
Antibody Diluent 4BR	1 x 6ml
Mouse IFN beta (Interferon beta 1) Lyophilized Recombinant Protein	2 vials
Plate Seals	1 unit
Sample Diluent NS (ab193972)	1 x 12ml
SimpleStep Pre-Coated 96-Well Microplate (ab206978)	1 unit
Stop Solution	1 x 12ml
TMB Development Solution	1 x 12ml

功能

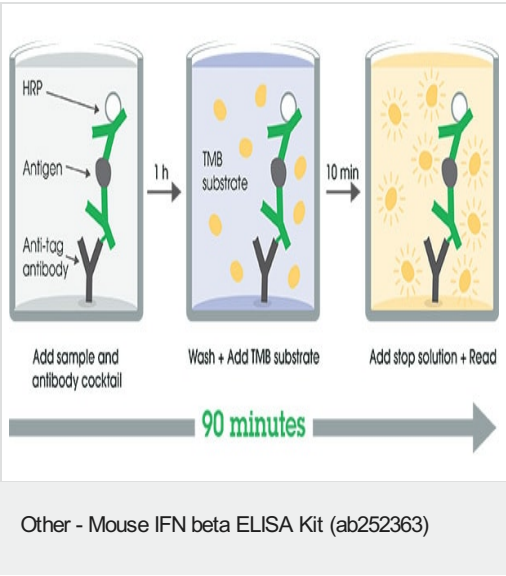
Has antiviral, antibacterial and anticancer activities.

序列相似性

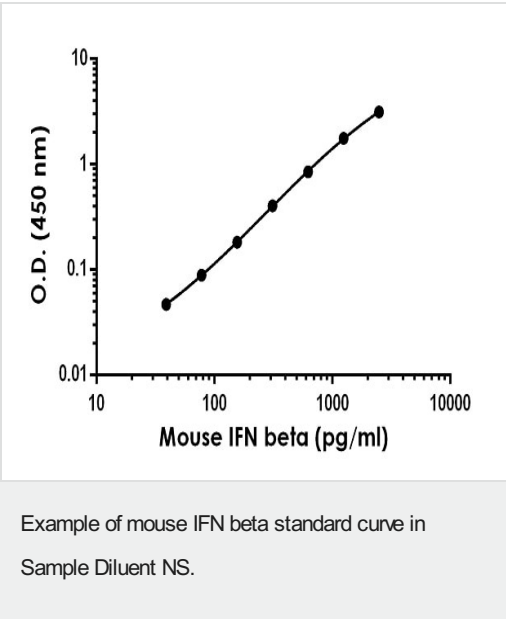
Belongs to the alpha/beta interferon family.

细胞定位

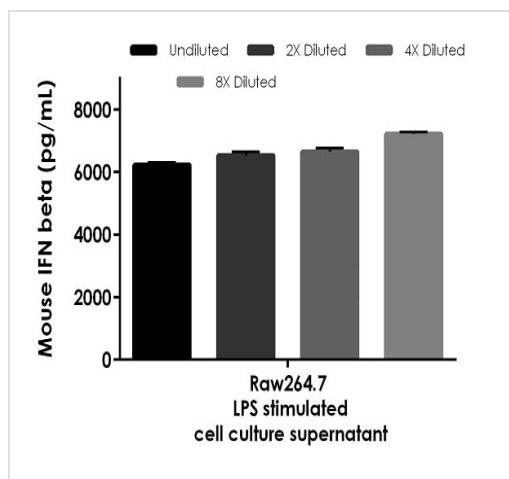
Secreted.



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.



The IFN beta 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.



Interpolated concentrations of native IFN beta in LPS stimulated Raw264.7 cell culture supernatant samples.

The concentrations of IFN beta were measured in duplicates, interpolated from the IFN beta standard curves and corrected for sample dilution. Undiluted samples are as follows: Raw264.7 cell culture supernatant 25%, The interpolated dilution factor corrected values are plotted (mean \pm SD, $n=2$). The mean IFN beta concentration was determined to be 6,664 pg/mL in Raw264.7 cell culture supernatant. Mouse IFN beta was below the detection point in unstimulated Raw264.7 supernatant.

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 - Mouse IFN beta ELISA Kit (ab252363)

To learn more about the advantages of recombinant antibodies 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.cn/abpromise> or contact our technical team.

Terms and conditions

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