

Human Follistatin ELISA Kit ab192147

SimpleStep ELISA

★★★★★ [1 Abreviews](#) [1 References](#) [6 图像](#)

概述

产品名称 人Follistatin ELISA试剂盒

检测方法 Colorimetric

批次内				
样品	n	Mean	SD	CV%
Overall	8			6.3%

批次间				
样品	n	Mean	SD	CV%
Overall	3			2.5%

样品类型 Cell culture supernatant, Serum, EDTA Plasma

检测类型 Sandwich (quantitative)

灵敏度 28 pg/ml

范围 0.08 ng/ml - 5 ng/ml

特定样本回收率		
样品类型	平均%	范围
Serum	110	99% - 125%
Cell culture media	86	85% - 86%
EDTA Plasma	92	89% - 95%

检测时间 1h 30m

实验步骤 One step assay

种属反应性 与反应: Human

产品概述 Human Follistatin ELISA Kit (ab192147) is a single-wash 90 min sandwich ELISA designed for the quantitative measurement of Follistatin protein in cell culture supernatant, edta plasma, and

serum. It uses our proprietary SimpleStep ELISA® technology. Quantitate Human Follistatin 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.

说明

Follistatin is a single-chain polypeptide consisting of about 344 amino acids. The follistatin protein is encoded by a relatively small gene size of 6 kb and 6 exons with alternative splice variants and post translational modifications to give a variety of molecular weights between 31 – 39 kDa. Follistatin forms a high affinity complex with activin sequestering it from its receptor thus having an inhibitory effect. Follistatin acts to promote muscle growth in mammals by acting as an antagonist to the protein myostatin. Myostatin acts as an inhibitor to muscle growth.

平台

Microplate

性能

存放说明

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

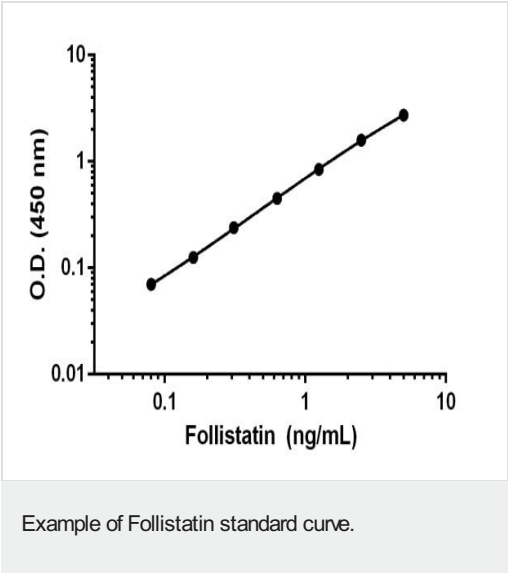
组件	1 x 96 tests
10X Human Follistatin Capture Antibody	1 x 600µl
10X Human Follistatin Detector Antibody	1 x 600µl
10X Wash Buffer PT (ab206977)	1 x 20ml
Antibody Diluent 5BI	1 x 6ml
Human Follistatin Lyophilized Recombinant Protein	2 vials
Plate Seals	1 unit
Sample Diluent 25BP	1 x 20ml
Sample Diluent NS (ab193972)	1 x 50ml
SimpleStep Pre-Coated 96-Well Microplate (ab206978)	1 unit
Stop Solution	1 x 12ml

组件	1 x 96 tests
TMB Development Solution	1 x 12ml

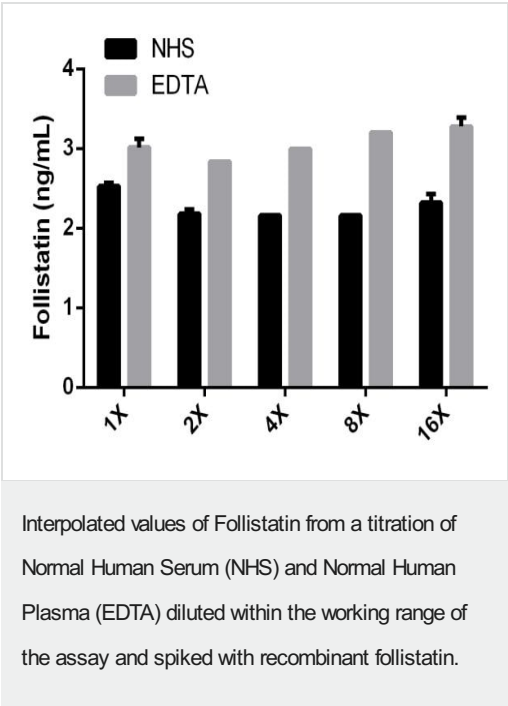
细胞定位

Secreted

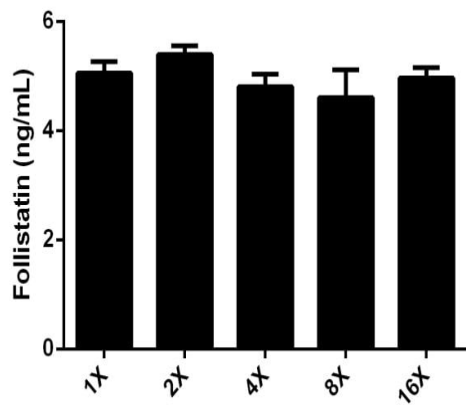
图片



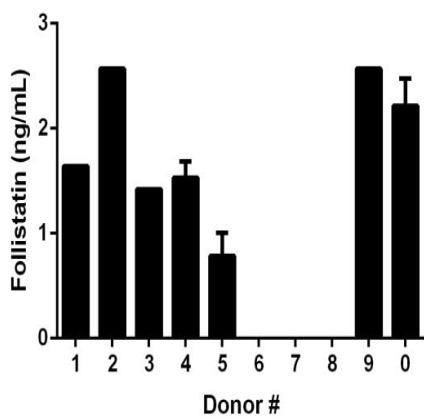
Background-subtracted data values (mean +/- SD) are graphed.



(mean +/- SD, n=2).

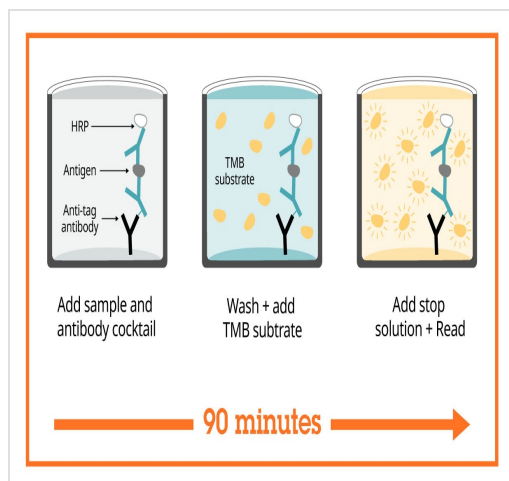


Interpolated values of Follistatin from HepG2 cell culture supernatants diluted within the working range of the assay (+/- SD, n=2).



Quantitation of Follistatin expression in serum samples diluted to within working range from individual donors.

Interpolated values of Follistatin are plotted on the bar graph. No signal was detected in Donors 6, 7 or 8.



Sandwich ELISA - Human Follistatin ELISA Kit
(ab192147)

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.

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 Follistatin ELISA Kit
(ab192147)

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

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Promise 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