

Human FGF19 ELISA Kit ab230943

重组 SimpleStep ELISA

8 图像

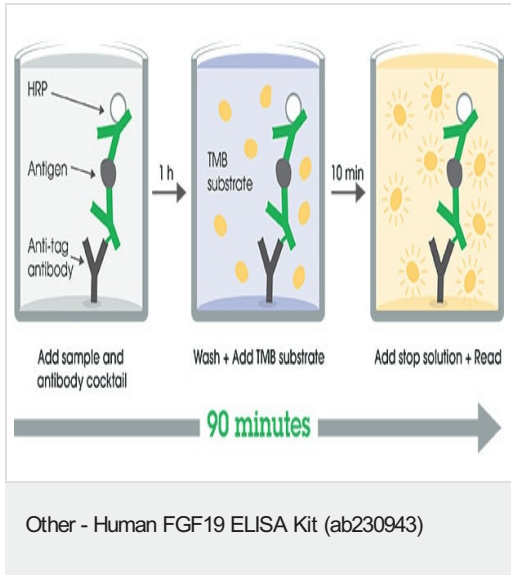
概述

产品名称	人FGF19 ELISA试剂盒				
检测方法	Colorimetric				
精确度	批次内				
样品		n	Mean	SD	CV%
Supernatant		5			3.8%
					批次间
样品		n	Mean	SD	CV%
Supernatant		3			6.6%
样品类型	Cell culture supernatant, Serum, Hep Plasma, EDTA Plasma, Cit plasma				
检测类型	Sandwich (quantitative)				
灵敏度	1.4 pg/ml				
范围	3.91 pg/ml - 250 pg/ml				
回收率	特定样本回收率				
样品类型		平均%		范围	
Serum		112		106% - 117%	
Cell culture media		108		107% - 110%	
Hep Plasma		108		101% - 116%	
EDTA Plasma				94% - 99%	
Cit plasma		100		97% - 105%	
检测时间	1h 30m				
实验步骤	One step assay				

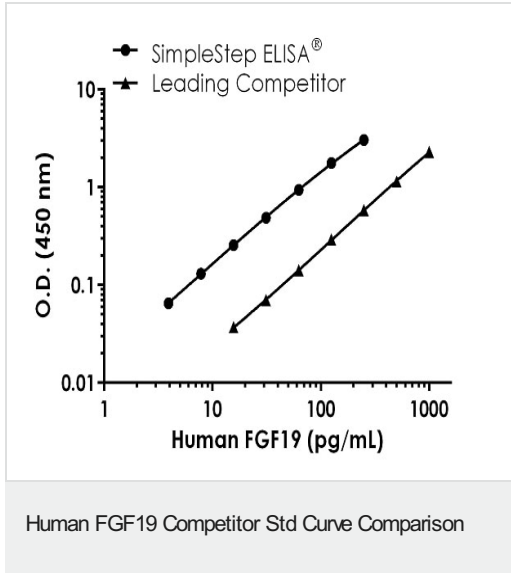
种属反应性	<p>与反应: Human</p> <p>不与反应: Cow</p>
产品概述	<p>Human FGF19 ELISA Kit (ab230943) is a single-wash 90 min sandwich ELISA designed for the quantitative measurement of FGF19 protein in cit plasma, edta plasma, hep plasma, and serum. It uses our proprietary SimpleStep ELISA® technology. Quantitate Human FGF19 with 1.4 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"> - 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 <p>A 384-well SimpleStep ELISA® microplate (ab203359) is available to use as an alternative to the 96-well microplate provided with SimpleStep ELISA® kits.</p>
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 Human FGF19 Capture Antibody	1 x 600µl
10X Human FGF19 Detector Antibody	1 x 600µl
10X Wash Buffer PT (ab206977)	1 x 20ml
Antibody Diluent 4BI	1 x 6ml
Human FGF19 Lyophilized Recombinant Protein	2 vials
Plate Seals	1 unit
Sample Diluent NS (ab193972)	1 x 50ml
SimpleStep Pre-Coated 96-Well Microplate (ab206978)	1 unit
Stop Solution	1 x 12ml
TMB Development Solution	1 x 12ml
功能	Involved in the suppression of bile acid biosynthesis through down-regulation of CYP7A1

	expression, following positive regulation of the JNK and ERK1/2 cascades. Stimulates glucose uptake in adipocytes. Activity requires the presence of KLB.
组织特异性	Expressed in fetal brain, cartilage, retina, and adult gall bladder.
序列相似性	Belongs to the heparin-binding growth factors family.
细胞定位	Secreted.

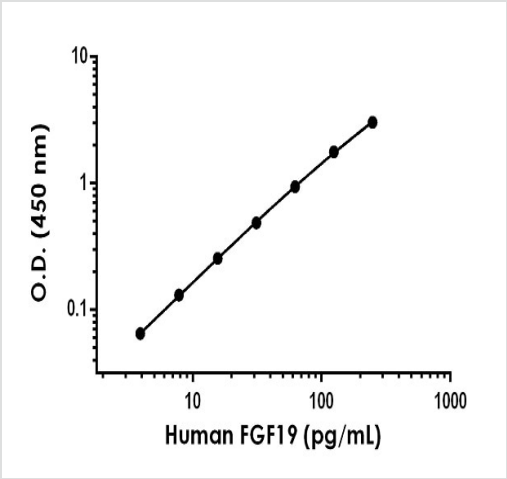
图片



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.

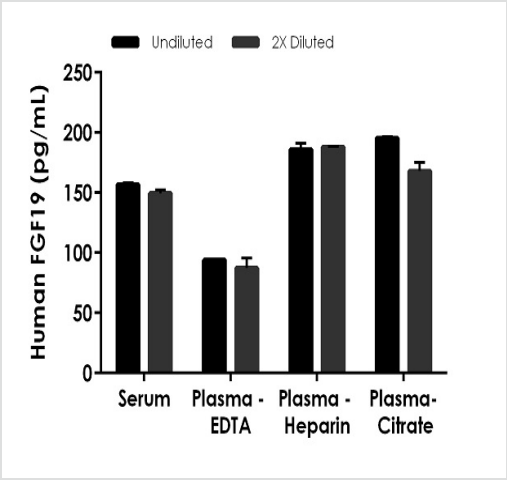


Standard curve comparison between human FGF19 SimpleStep ELISA® kit and traditional ELISA kit from leading competitor. SimpleStep ELISA kit shows increased sensitivity.



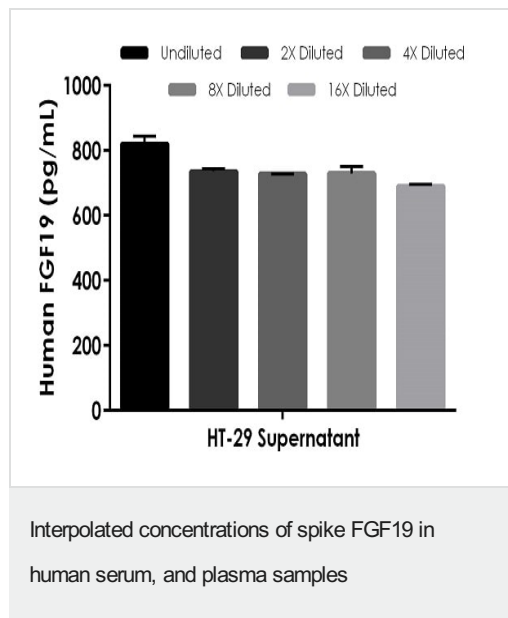
Background-subtracted data values (mean +/- SD) are graphed.

Example of human FGF19 standard curve in Sample Diluent NS

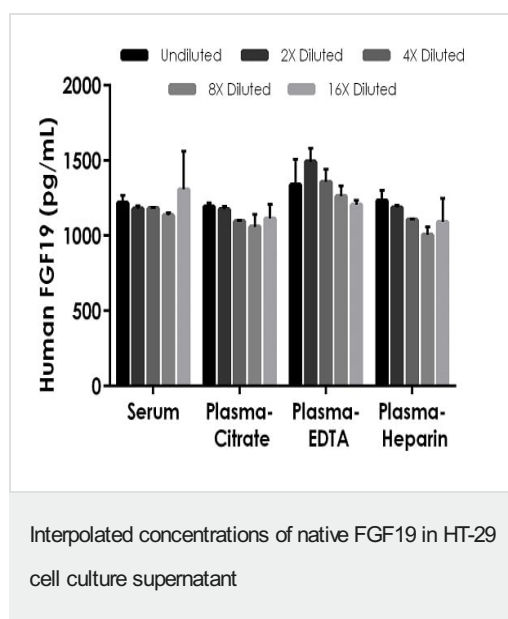


The concentrations of FGF19 were measured in duplicates, interpolated from the FGF19 standard curves and corrected for sample dilution. Undiluted samples are as follows: serum 12.5%, plasma (EDTA) 12.5%, plasma (heparin) 12.5% and plasma (citrate) 12.5%. The interpolated dilution factor corrected values are plotted (mean +/- SD, n=2). The mean FGF19 concentration was determined to be 153 pg/mL in serum , 91 pg/mL in plasma (EDTA), 187 pg/mL in plasma (heparin), and 182 pg/mL in plasma (citrate).

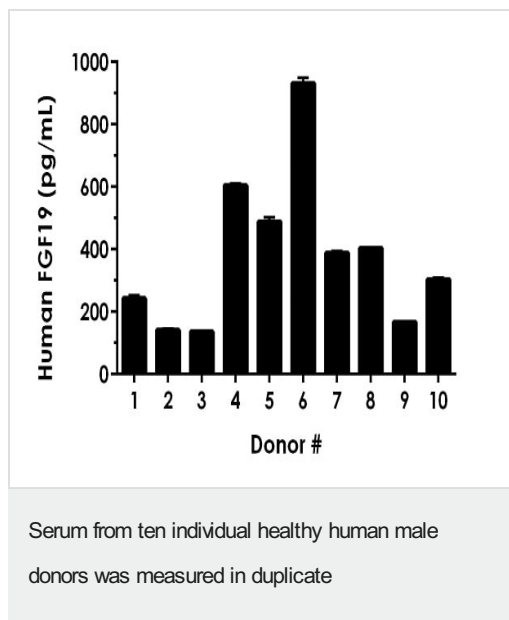
Interpolated concentrations of native FGF19 in human serum, and plasma samples



The concentrations of FGF19 were measured in duplicates, interpolated from the FGF19 standard curves and corrected for sample dilution. Undiluted samples are as follows: serum 12.5%, plasma (citrate) 12.5%, plasma (EDTA) 12.5% and plasma (heparin) 12.5%. The interpolated dilution factor corrected values are plotted (mean \pm SD, n=2).



The concentrations of FGF19 were measured in duplicates, interpolated from the FGF19 standard curves and corrected for sample dilution. Undiluted samples are as follows: HT-29 cell culture supernatant 25%. The interpolated dilution factor corrected values are plotted (mean \pm SD, n=2). The mean FGF19 concentration was determined to be 741 pg/mL in HT-29 cell culture supernatant.



Interpolated dilution factor corrected values are plotted (mean \pm SD, $n=2$). The mean FGF19 concentration was determined to be 381 pg/mL with a range of 136 – 944 pg/mL.

Powered by
recombinant antibodies

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

Sandwich ELISA - Human FGF19 ELISA Kit
(ab230943)

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