

Human aFGF ELISA Kit (acid FGF) ab193691

[1 References](#) [1 图像](#)

概述

产品名称 人aFGF ELISA试剂盒(acid FGF)

检测方法 Colorimetric

精确度 批次内

样品	n	Mean	SD	CV%
Overall				< 10%

批次间

样品	n	Mean	SD	CV%
Overall				< 12%

样品类型 Cell culture supernatant, Serum, Plasma

检测类型 Sandwich (quantitative)

灵敏度 12 pg/ml

范围 12 pg/ml - 4000 pg/ml

回收率 特定样本回收率

样品类型	平均%	范围
Serum	122.5	105% - 131%
Plasma	91.51	78% - 103%
Cell culture media	76.17	66% - 87%

实验步骤 Multiple steps standard assay

种属反应性 与反应: Human

产品概述 Abcam's aFGF (acid FGF) Human ELISA Kit (ab193691) is an in vitro enzyme-linked immunosorbent assay for the quantitative measurement of Human aFGF in serum, plasma and cell culture supernatants.

This assay employs an antibody specific for human aFGF coated on a 96-well plate. Standards and samples are pipetted into the wells and aFGF present in a sample is bound to the wells by

the immobilized antibody. The wells are washed and biotinylated anti-human aFGF antibody is added. After washing away any unbound biotinylated antibody, an HRP-conjugated streptavidin is pipetted to the wells. After incubation, the wells are again washed, followed by the addition of a TMB substrate solution to the wells. Color will develop in proportion to the amount of aFGF bound in each well. Addition of the Stop Solution will change the color from blue to yellow, and the intensity of the color is measured at 450 nm.

平台 Pre-coated microplate (12 x 8 well strips)

性能

存放说明 Store at -20°C. Please refer to protocols.

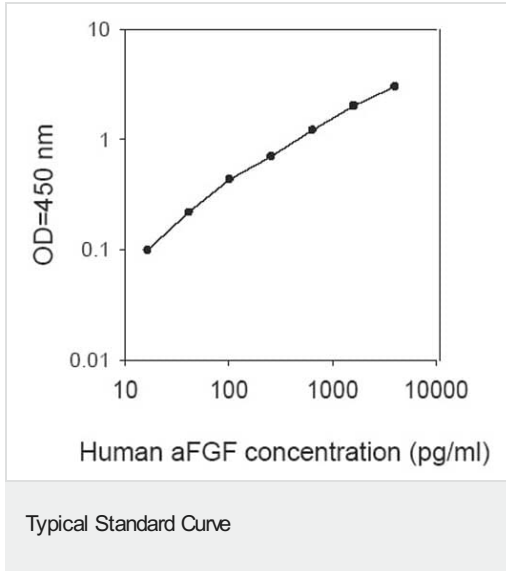
组件	1 x 96 tests
20X Wash Buffer	1 x 25ml
5X Assay Diluent	1 x 15ml
500X HRP-Streptavidin Concentrate	1 x 200µl
Biotinylated Human aFGF detection antibody	2 vials
Human aFGF standards (lyophilized)	2 vials
Pre-coated Human aFGF Microplate (12 strips x 8 wells)	1 unit
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

功能 The heparin-binding fibroblast growth factors play important roles in the regulation of cell survival, cell division, angiogenesis, cell differentiation and cell migration. They are potent mitogens in vitro.

序列相似性 Belongs to the heparin-binding growth factors family.

细胞定位 Secreted. Cytoplasm. Cytoplasm > cell cortex. Lacks a cleavable signal sequence. Within the cytoplasm, it is transported to the cell membrane and then secreted by a non-classical pathway that requires Cu(2+) ions and S100A13. Secreted in a complex with SYT1.

图片



Standard Curve using ab193691 aFGF (acid FGF) Human ELISA Kit.

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