

Human αFGF ELISA Kit (acid FGF) ab193691

1 References 1 Image

Overview

Product name Human αFGF ELISA Kit (acid FGF)

Detection method Colorimetric

Precision Intra-assay

Sample	n	Mean	SD	CV%
Overall				< 10%

Inter-assay

Sample	n	Mean	SD	CV%
Overall				< 12%

Sample type Cell culture supernatant, Serum, Plasma

Assay type Sandwich (quantitative)

Sensitivity 12 pg/ml

Range 12 pg/ml - 4000 pg/ml

Recovery Sample specific recovery

Sample type	Average %	Range
Serum	122.5	105% - 131%
Plasma	91.51	78% - 103%
Cell culture media	76.17	66% - 87%

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview Abcam’s αFGF (acid FGF) Human ELISA Kit (ab193691) is an in vitro enzyme-linked immunosorbent assay for the quantitative measurement of Human αFGF in serum, plasma and cell culture supernatants.

This assay employs an antibody specific for human αFGF coated on a 96-well plate. Standards and samples are pipetted into the wells and αFGF 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.

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

Properties

Storage instructions Store at -20°C. Please refer to protocols.

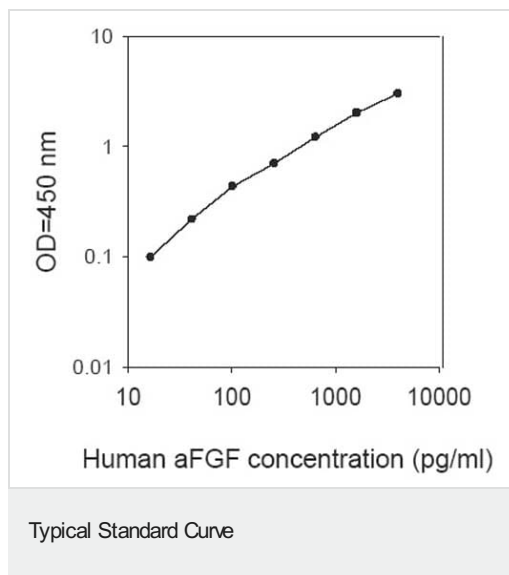
Components	1 x 96 tests
20X Wash Buffer	1 x 25ml
5X Assay Diluent	1 x 15ml
500X HRP-Streptavidin Concentrate	1 x 200µl
Biotinylated Human aFGF detection antibody	2 vials
Human aFGF standards (lyophilized)	2 vials
Pre-coated Human aFGF Microplate (12 strips x 8 wells)	1 unit
Stop Solution	1 x 8ml
TMB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

Function 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.

Sequence similarities Belongs to the heparin-binding growth factors family.

Cellular localization 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.

Images



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.com/abpromise> or contact our technical team.

Terms and conditions

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