

Product datasheet

Human FGF4 ELISA Kit ab100517

[1 Image](#)

Overview

Product name	Human FGF4 ELISA Kit
Detection method	Colorimetric
Sample type	Cell culture extracts, Tissue Extracts
Assay type	Sandwich (quantitative)
Sensitivity	< 50 pg/ml
Range	24.69 pg/ml - 18000 pg/ml
Recovery	92 %

Sample specific recovery

Sample type	Average %	Range
Cell culture extracts	93.45	94% - 103%
Tissue Extracts	91.96	82% - 102%

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview Abcam's FGF4 Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Human FGF4 in cell lysates and tissue lysates.

This assay employs an antibody specific for Human FGF4 coated on a 96-well plate. Standards and samples are pipetted into the wells and FGF4 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human FGF4 antibody is added. After washing away unbound biotinylated antibody, HRP-conjugated streptavidin is pipetted to the wells. The wells are again washed, a TMB substrate solution is added to the wells and color develops in proportion to the amount of FGF4 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

Tested applications **Suitable for:** Sandwich ELISA

Platform Microplate

Properties

Storage instructions

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

Components	1 x 96 tests
120X HRP-Streptavidin Concentrate	1 x 200µl
20X Wash Buffer Concentrate	1 x 25ml
2x Cell lysate buffer	1 x 5ml
5X Assay Diluent	1 x 15ml
5X Sample Diluent Buffer	1 x 10ml
Biotinylated anti-Human FGF4	2 vials
FGF4 Microplate (12 strips x 8 wells)	1 unit
Recombinant Human FGF4 Standard	2 vials
Stop Solution	1 x 8ml
TMB Substrate Reagent	1 x 12ml

Relevance

FGF4 is a heparin binding growth factor that is a member of the FGF family. Proteins of this family play a central role during prenatal development and postnatal growth and regeneration of a variety of tissues, by promoting cellular proliferation and differentiation. FGF4 signals through the FGFR 1c, 2c, 3c, and 4.

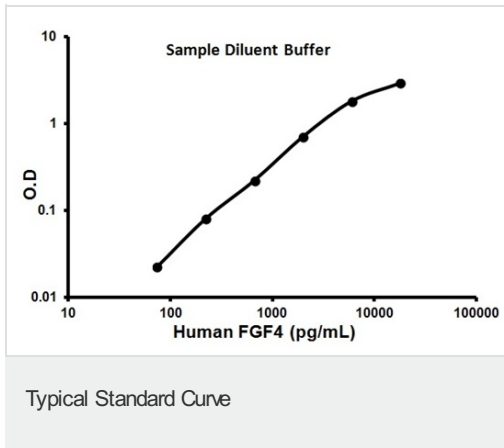
Applications

Our [Abpromise guarantee](#) covers the use of **ab100517** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

Application	Abreviews	Notes
Sandwich ELISA		Use at an assay dependent concentration.

Images



Representative standard curve using
ab100517

Please note: All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

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 <http://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