

Human FGF basic ELISA Kit (FGF2) αb99980

1 References 3 Images

Overview

Product name	Human FGF basic ELISA Kit (FGF2)
Detection method	Colorimetric
Sample type	Cell culture extracts, Tissue Extracts
Assay type	Sandwich (quantitative)
Sensitivity	< 50 pg/ml
Range	102.4 pg/ml - 10000 pg/ml
Recovery	> 95 %

Sample specific recovery

Sample type	Average %	Range
Cell culture extracts	97.65	87% - 108%
Tissue Extracts	95.87	86% - 107%

Assay duration	Multiple steps standard assay
Species reactivity	Reacts with: Human
Product overview	<p>Abcam's FGF basic (FGF2) Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an <i>in vitro</i> enzyme-linked immunosorbent assay for the quantitative measurement of Human FGF basic in cell lysates and tissue lysates.</p> <p>This assay employs an antibody specific for Human FGF basic coated on a 96-well plate. Standards and samples are pipetted into the wells and FGF basic present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human FGF basic 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 FGF basic bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.</p>

Notes	<p>Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances.</p> <p>It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.</p>
-------	--

Platform

Microplate

Properties**Storage instructions**

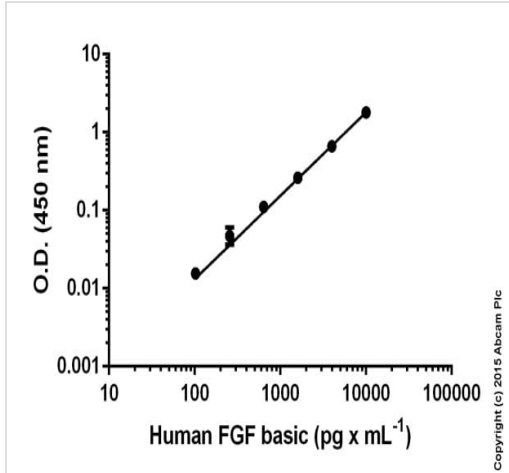
Store at -20°C. Please refer to protocols.

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

Relevance

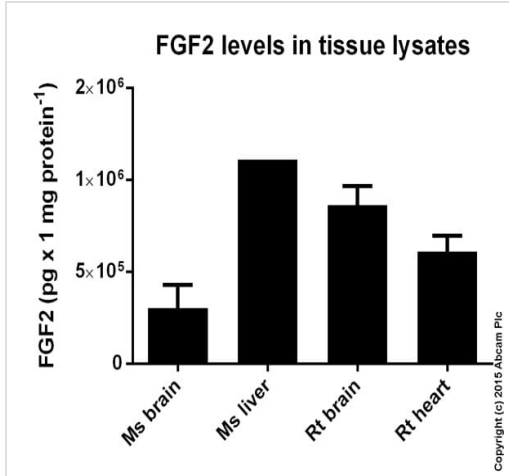
FGF basic (FGF2) is a member of the fibroblast growth factor (FGF) family. FGF family members bind heparin and possess broad mitogenic and angiogenic activities. FGF2 is involved in diverse biological processes, such as limb and nervous system development, wound healing, and tumor growth. FGF2 mRNA contains multiple polyadenylation sites, and is alternatively translated from AUG and non-AUG (CUG) initiation codons resulting in five different isoforms with distinct properties. The CUG-initiated isoforms are localized in the nucleus and are responsible for the intracrine effect, whereas, the AUG-initiated form is mostly cytosolic and is responsible for the paracrine and autocrine effects of this FGF.

Images



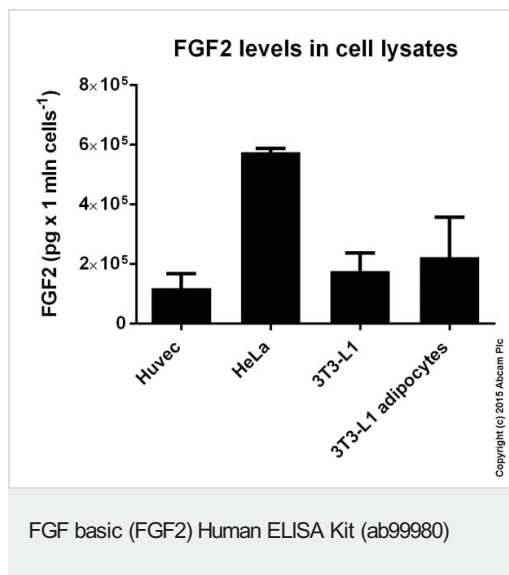
Standard curve: mean of duplicates (+/- SD) with background reads subtracted

FGF basic (FGF2) Human ELISA Kit (ab99980)



FGF2 measured in tissue lysates showing quantity (pg) per 1 mg of extracted protein. Protein concentration for samples varied from 7 mg/mL to 17 mg/mL. Samples were diluted 200-2000 fold.

FGF basic (FGF2) Human ELISA Kit (ab99980)



FGF2 measured in cell culture lysates showing quantity (pg) per 1 mln cells. Samples with the concentration of 1e7 cells/mL were used. Samples were diluted 200-2000 fold.

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