

Product datasheet

Human FGF23 ELISA Kit ab267652

1 Image

Overview

**Product name** Human FGF23 ELISA Kit

**Detection method** Colorimetric

**Precision**

Intra-assay

Sample	n	Mean	SD	CV%
Overall				< 10%

Inter-assay

Sample	n	Mean	SD	CV%
Overall				< 12%

**Sample type** Serum, Plasma, Cell culture media

**Assay type** Sandwich (quantitative)

**Sensitivity** 0.3 ng/ml

**Range** 0.307 ng/ml - 75 ng/ml

**Recovery**

Sample specific recovery

Sample type	Average %	Range
Serum	135	122% - 145%
Plasma	144	138% - 147%
Cell culture media	143	133% - 149%

**Assay duration** Multiple steps standard assay

**Species reactivity** **Reacts with:** Human

**Product overview** Human FGF23 ELISA Kit is designed for the quantitative determination of FGF23 in cell culture supernatants, plasma and serum samples.

**ΔNote:** Human FGF-23 concentration is low in normal serum/plasma, and may not be detectable in this assay.

This assay employs an antibody specific for human FGF23 coated on a 96-well plate. Standards and samples are pipetted into the wells and FGF23 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-human FGF23 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 FGF23 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

**Platform** Microplate

## Properties

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

Components	1 x 96 tests
20X Wash Buffer	1 x 25ml
350X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent B	1 x 15ml
Anti-Human FGF23 coated Microplate (12 x 8 wells)	1 unit
Assay Diluent C	1 x 30ml
Biotinylated Anti-Human FGF23 Detection Antibody	2 vials
Human FGF23 Standard (Lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB Substrate Solution	1 x 12ml

**Function** Regulator of phosphate homeostasis. Inhibits renal tubular phosphate transport by reducing SLC34A1 levels. Upregulates EGR1 expression in the presence of KL (By similarity). Acts directly on the parathyroid to decrease PTH secretion (By similarity). Regulator of vitamin-D metabolism. Negatively regulates osteoblast differentiation and matrix mineralization.

**Tissue specificity** Expressed in osteogenic cells particularly during phases of active bone remodeling. In adult trabecular bone, expressed in osteocytes and flattened bone-lining cells (inactive osteoblasts).

**Involvement in disease** Defects in FGF23 are the cause of autosomal dominant hypophosphataemic rickets (ADHR) [MIM:193100]. ADHR is characterized by low serum phosphorus concentrations, rickets, osteomalacia, leg deformities, short stature, bone pain and dental abscesses. Defects in FGF23 are a cause of hyperphosphatemic familial tumoral calcinosis (HFTC) [MIM:211900]. HFTC is a severe autosomal recessive metabolic disorder that manifests with hyperphosphatemia and massive calcium deposits in the skin and subcutaneous tissues.

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

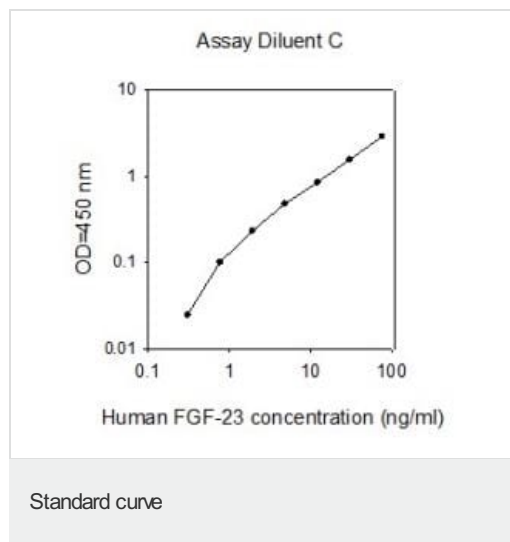
**Post-translational modifications** Following secretion this protein is inactivated by cleavage into a N-terminal fragment and a C-terminal fragment. The processing is effected by proprotein convertases. O-glycosylated by GALT3. Glycosylation is necessary for secretion; it blocks processing by

proprotein convertases when the O-glycan is alpha 2,6-sialylated. Competition between proprotein convertase cleavage and block of cleavage by O-glycosylation determines the level of secreted active FGF23.

#### Cellular localization

Secreted. Secretion is dependent on O-glycosylation.

#### Images



Human FGF23 ELISA kit (ab267652) Standard curve.

Data provided for demonstration purposes only. A new standard curve must be generated for each assay performed.

**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