

### Human GDF-15 ELISA Kit ab155432

★★★★★ [2 Abreviews](#) [6 References](#) [3 Images](#)

#### Overview

**Product name** Human GDF-15 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** Cell culture supernatant, Serum, Plasma

**Assay type** Sandwich (quantitative)

**Sensitivity** < 2 pg/ml

**Range** 1.1 pg/ml - 800 pg/ml

**Recovery** Sample specific recovery

Sample type	Average %	Range
Serum	104.2	92% - 116%
Plasma	102.9	88% - 119%
Cell culture media	114.2	96% - 105%

**Assay duration** Multiple steps standard assay

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

**Predicted to work with:** Non human primates 

**Product overview** Abcam's GDF-15 Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay designed for the quantitative measurement of Human GDF-15 in serum, plasma and cell culture supernatants.

This assay employs an antibody specific for Human GDF-15 coated on a 96-well plate. Standards and samples are pipetted into the wells and GDF-15 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human GDF-15 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 GDF-15 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.

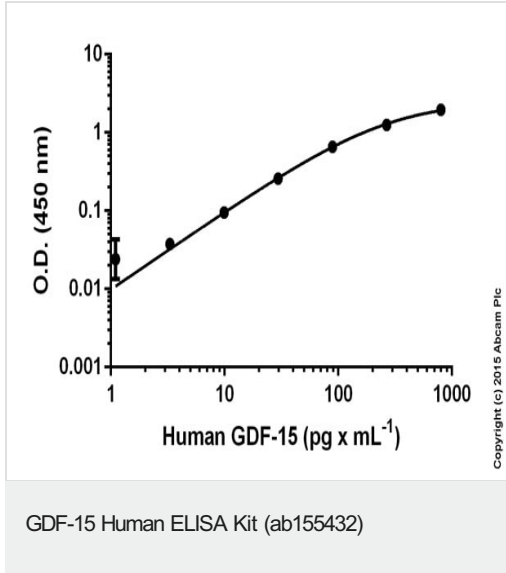
Components	1 x 96 tests
20X Wash Buffer	1 x 25ml
5X Assay Diluent	1 x 15ml
500X HRP-Streptavidin Concentrate	1 x 200µl
Biotinylated anti-Human GDF-15	2 vials
GDF-15 Microplate (12 x 8 wells)	1 unit
Recombinant Human GDF-15 Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

**Tissue specificity** Highly expressed in placenta, with lower levels in prostate and colon and some expression in kidney.

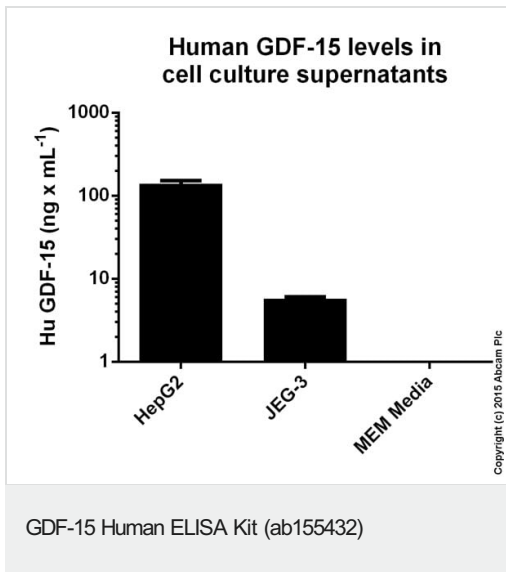
**Sequence similarities** Belongs to the TGF-beta family.

**Cellular localization** Secreted.

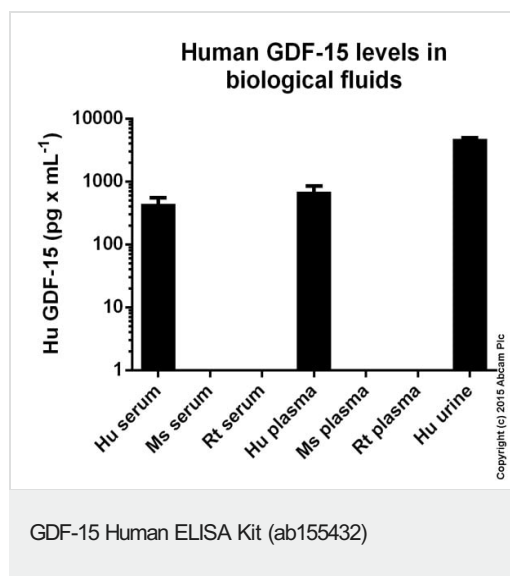
## Images



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



GDF-15 measured in cell culture supernatants showing quantity (ng) per mL of tested sample. Samples were diluted 20-540 fold.



GDF-15 measured in biological fluids showing quantity (pg) per mL of tested sample. Samples were diluted 2-54 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