

Human IGF1 ELISA Kit ab100545

★★★★★ [2 Abreviews](#) [23 References](#) [画像数 3](#)

製品の概要

| | | | | | |
|---------|--|-------|-----|-------------|-------|
| 製品名 | Human IGF1 ELISA Kit | | | | |
| 検出方法 | Colorimetric | | | | |
| 再現性 | Intra-Assay (同時再現性) | | | | |
| | サンプル | N | 平均値 | SD | CV% |
| | Overall | | | | < 10% |
| | Inter-Assay (日差再現性) | | | | |
| | サンプル | N | 平均値 | SD | CV% |
| | Overall | | | | < 12% |
| サンプルの種類 | Cell culture supernatant, Serum, Plasma | | | | |
| アッセイタイプ | Sandwich (quantitative) | | | | |
| 検出感度 | < 0.2 ng/ml | | | | |
| 検出範囲 | 0.1 ng/ml - 30 ng/ml | | | | |
| 添加回収試験 | 97 % | | | | |
| | 特定サンプルでの回収試験 | | | | |
| | サンプルの種類 | 平均 % | | 測定範囲 | |
| | Cell culture supernatant | 102.5 | | 98% - 110% | |
| | Serum | 119.6 | | 105% - 138% | |
| | Plasma | 105.4 | | 81% - 121% | |
| ステップ | Multiple steps standard assay | | | | |
| 種交差性 | 交差種: Human | | | | |
| 製品の概要 | Abcam's IGF1 Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an <i>in vitro</i> enzyme linked immunosorbent assay for the quantitative measurement of Human IGF1 in serum, plasma, and cell culture supernatants. | | | | |

This ELISA kit can only detect the free form of human IGF-1. It does not detect the complex or bound forms such as IGFBP-bound IGF1.

This assay employs an antibody specific for Human IGF1 coated on a 96-well plate. Standards and samples are pipetted into the wells and IGF1 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human IGF1 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 IGF1 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

Get results in 90 minutes with Human IGF1 ELISA Kit ([ab211651](#)) from our SimpleStep ELISA® range.

特記事項

Optimization may be required with urine samples

試験プラットフォーム

Microplate

製品の特性

保存方法

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

| 内容 | 1 x 96 tests |
|---|--------------|
| 120X HRP-Streptavidin Concentrate | 1 x 200µl |
| 20X Wash Buffer | 1 x 25ml |
| Assay Diluent C | 2 x 30ml |
| Biotinylated anti-Human IGF-1 | 2 vials |
| IGF-1 Microplate (12 strips x 8 wells) | 1 unit |
| Recombinant Human IGF1 Standard (lyophilized) | 2 vials |
| Stop Solution | 1 x 8ml |
| TMB One-Step Substrate Reagent | 1 x 12ml |

機能

The insulin-like growth factors, isolated from plasma, are structurally and functionally related to insulin but have a much higher growth-promoting activity. May be a physiological regulator of [1-14C]-2-deoxy-D-glucose (2DG) transport and glycogen synthesis in osteoblasts. Stimulates glucose transport in rat bone-derived osteoblastic (PyMS) cells and is effective at much lower concentrations than insulin, not only regarding glycogen and DNA synthesis but also with regard to enhancing glucose uptake.

関連疾患

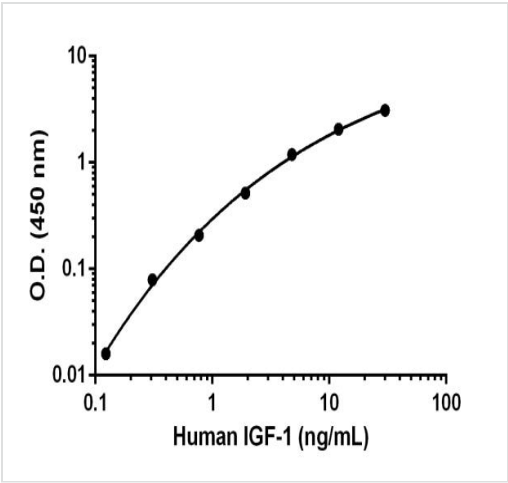
Defects in IGF1 are the cause of insulin-like growth factor I deficiency (IGF1 deficiency) [MIM:608747]. IGF1 deficiency is an autosomal recessive disorder characterized by growth retardation, sensorineural deafness and mental retardation.

配列類似性

Belongs to the insulin family.

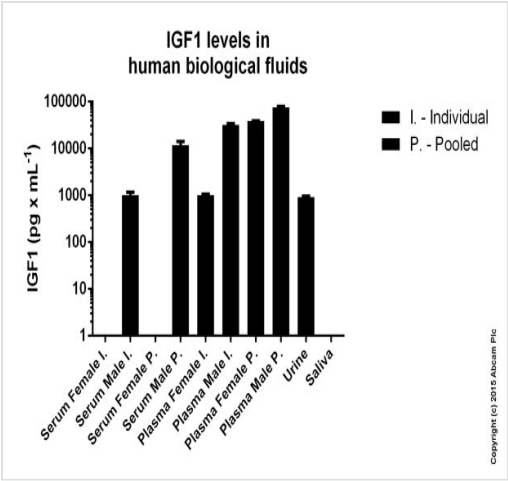
細胞内局在

Secreted.



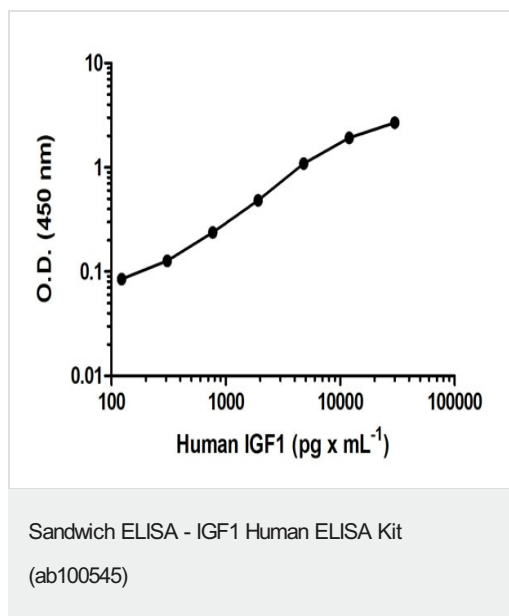
Typical Standard Curve

Representative standard curve using ab100545.



Sandwich ELISA - IGF1 Human ELISA Kit
(ab100545)

IGF1 measured in biological fluids showing quantity (pg) per mL of tested sample. Samples were diluted 1-64 fold.



Human IGF1 standard curve: mean of duplicates (+/- SD)

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.co.jp/abpromise> or contact our technical team.

Terms and conditions

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