

Mouse IGF1 ELISA Kit ab100695

★★★★★ [1 Abreviews](#) [31 References](#) [画像数 4](#)

製品の概要

製品名	Mouse IGF1 ELISA Kit
検出方法	Colorimetric
サンプルの種類	Cell culture supernatant, Serum, Plasma, Tissue Lysate
アッセイタイプ	Sandwich (quantitative)
検出感度	< 4 pg/ml
検出範囲	2.74 pg/ml - 2000 pg/ml
添加回収試験	98 %

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	86.4	72% - 99%
Serum	101.2	85% - 117%
Plasma	106.6	85% - 114%

ステップ Multiple steps standard assay

種交差性 交差種: Mouse

製品の概要

Abcam's IGF1 Mouse ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of mouse IGF1 in serum, plasma and cell culture supernatants.

This assay employs an antibody specific for mouse 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-mouse 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.

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
200X HRP-Streptavidin Concentrate	1 x 200µl
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
Assay Diluent A	1 x 30ml
Biotinylated anti-Mouse IGF1	2 vials
IGF1 Microplate (12 x 8 wells)	1 unit
Recombinant Mouse 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.

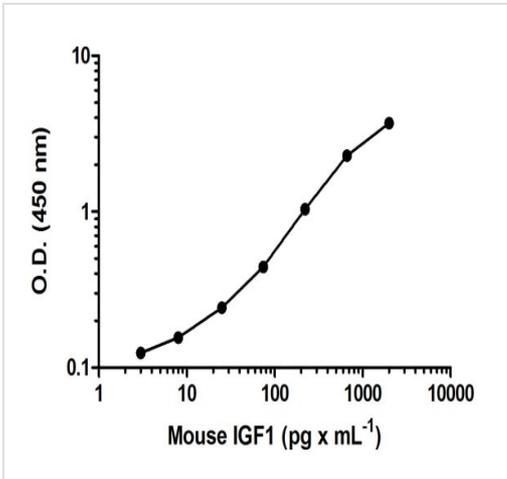
配列類似性

Belongs to the insulin family.

細胞内局在

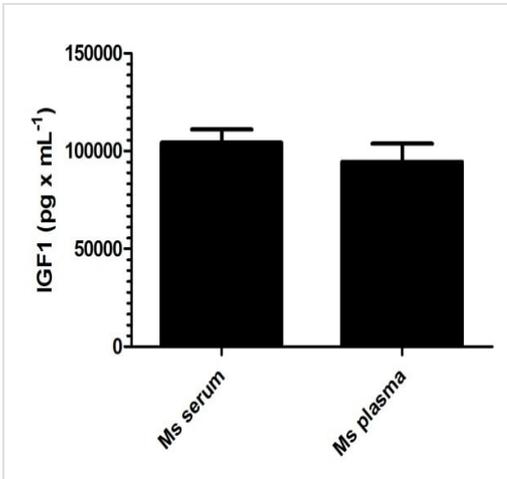
Secreted.

画像



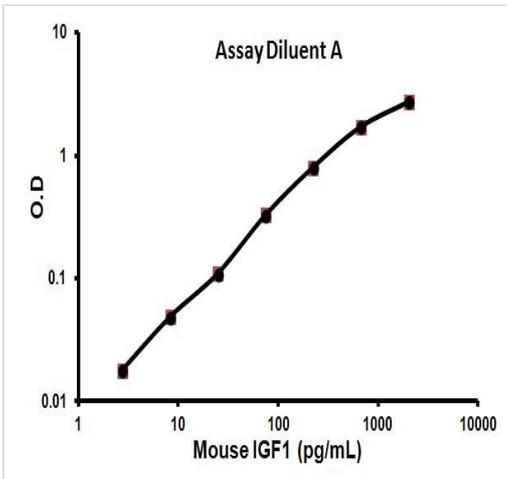
Mouse IGF1 standard curve: mean of duplicates (+/- SD), performed in Diluent A

Sandwich ELISA - IGF1 Mouse ELISA Kit
(ab100695)



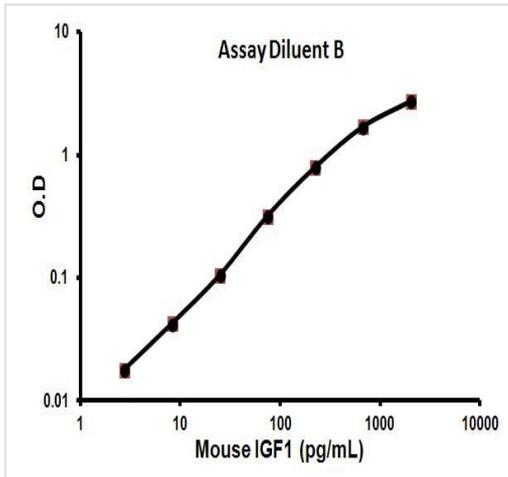
Mouse IGF1 measured in biological fluids showing quantity (pg) per mL of tested sample. Samples diluted 500-4500 fold.

Sandwich ELISA - IGF1 Mouse ELISA Kit
(ab100695)



Representative Standard Curve using ab100695

Typical Standard Curve



Representative Standard Curve using ab100695

Typical Standard Curve

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