

Mouse IGF2 ELISA Kit ab100696

[6 References](#) [1 Image](#)

Overview

| | |
|---------------------------|---|
| Product name | Mouse IGF2 ELISA Kit |
| Detection method | Colorimetric |
| Sample type | Cell culture supernatant |
| Assay type | Sandwich (quantitative) |
| Sensitivity | < 1.5 ng/ml |
| Range | 2.3 ng/ml - 150 ng/ml |
| Assay duration | Multiple steps standard assay |
| Species reactivity | Reacts with: Mouse |
| Product overview | <p>Abcam's IGF2 Mouse ELISA (Enzyme-Linked Immunosorbent Assay) kit is an <i>in vitro</i> enzyme-linked immunosorbent assay for the quantitative measurement of mouse IGF2 in cell culture supernatants.</p> <p>This assay employs an antibody specific for mouse IGF2 coated on a 96-well plate. Standards and samples are pipetted into the wells and IGF2 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-mouse IGF2 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 IGF2 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.</p> <p>Please note that ab100696 has only been validated with cell culture supernatant samples and cannot be used with plasma samples. It is also not validated for serum samples.</p> |
| Platform | Microplate |

Properties

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

| Components | 1 x 96 tests |
|-------------------------------------|--------------|
| 1,000X HRP-Streptavidin Concentrate | 1 x 200µl |
| 20X Wash Buffer | 1 x 25ml |

| Components | 1 x 96 tests |
|---|--------------|
| 5X Assay Diluent | 1 x 15ml |
| Biotinylated anti-Mouse IGF2 | 2 vials |
| IGF2 Microplate (12 x 8 wells) | 1 unit |
| Recombinant Mouse IGF2 Standard (lyophilized) | 2 vials |
| Stop Solution | 1 x 8ml |
| TMB One-Step Substrate Reagent | 1 x 12ml |

Function

The insulin-like growth factors possess growth-promoting activity. In vitro, they are potent mitogens for cultured cells. IGF-II is influenced by placental lactogen and may play a role in fetal development.

Preptin undergoes glucose-mediated co-secretion with insulin, and acts as physiological amplifier of glucose-mediated insulin secretion. Exhibits osteogenic properties by increasing osteoblast mitogenic activity through phosphoactivation of MAPK1 and MAPK3.

Involvement in disease

Epigenetic changes of DNA hypomethylation in IGF2 are a cause of Silver-Russell syndrome (SIRS) [MIM:180860]. SIRS is a clinically heterogeneous condition characterized by severe intrauterine growth retardation, poor postnatal growth, craniofacial features such as a triangular shaped face and a broad forehead, body asymmetry, and a variety of minor malformations.

Sequence similarities

Belongs to the insulin family.

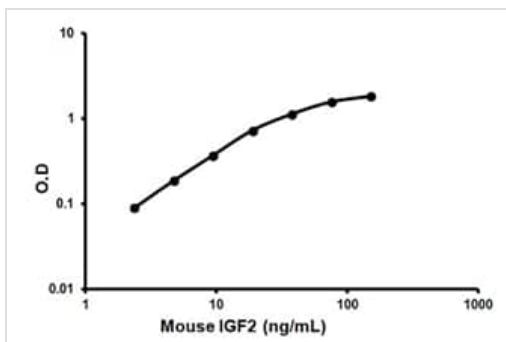
Post-translational modifications

O-glycosylated with a core 1 or possibly core 8 glycan.

Cellular localization

Secreted.

Images



Typical standard curve for ab100696

Standard Curve for Mouse IGF2

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