

Mouse IGF2 ELISA Kit ab100696

[6 References](#) [画像数 1](#)

製品の概要

製品名	Mouse IGF2 ELISA Kit
検出方法	Colorimetric
サンプルの種類	Cell culture supernatant
アッセイタイプ	Sandwich (quantitative)
検出感度	< 1.5 ng/ml
検出範囲	2.3 ng/ml - 150 ng/ml
ステップ	Multiple steps standard assay
種交差性	交差種: Mouse
製品の概要	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.

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.

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.

試験プラットフォーム Microplate

製品の特性

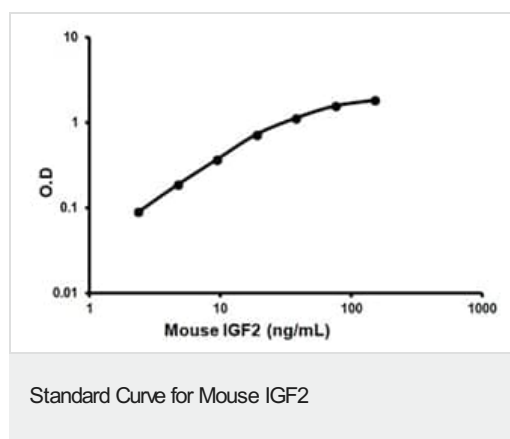
保存方法 Store at -20°C. Please refer to protocols.

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

内容	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

機能	<p>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.</p> <p>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.</p>
関連疾患	<p>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.</p>
配列類似性	<p>Belongs to the insulin family.</p>
翻訳後修飾	<p>O-glycosylated with a core 1 or possibly core 8 glycan.</p>
細胞内局在	<p>Secreted.</p>

画像



Typical standard curve for ab100696

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