

Mouse Insulin like Growth Factor 1 ELISA kit (IGF1) ab108874

5 References [画像数 1](#)

製品の概要

製品名 Mouse Insulin like Growth Factor 1 ELISA kit (IGF1)

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				3.2%

Inter-Assay (日差再現性)

サンプル	N	平均値	SD	CV%
Overall				9.8%

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 = 0.062 ng/ml

検出範囲 0.3 ng/ml - 5 ng/ml

添加回収試験 96 %

全工程の試験時間 5h 00m

ステップ Multiple steps standard assay

種交差性 交差種: Mouse

製品の概要 Abcam's Insulin like Growth Factor 1 (IGF1) Mouse *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of Insulin like Growth Factor 1 levels in plasma, serum and cell culture supernatants.

An IGF1 specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently an IGF1 specific biotinylated detection antibody is added and then followed by washing with wash buffer. Streptavidin-Peroxidase Conjugate is added and unbound conjugates are washed away with wash buffer. TMB is then used to visualize Streptavidin-Peroxidase enzymatic reaction. TMB is catalyzed by Streptavidin-Peroxidase to produce a blue color product that changes into yellow after adding acidic stop solution. The density of yellow coloration is directly proportional to the amount of IGF1 captured in plate.

The entire kit may be stored at -20°C for long term storage before reconstitution - Avoid repeated freeze-thaw cycles.

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
100X Streptavidin-Peroxidase Conjugate	1 x 80µl
10X Diluent N Concentrate	1 x 30ml
1X Standard Diluent	1 x 2ml
20X Wash Buffer Concentrate	2 x 30ml
40X Biotinylated Mouse Insulin like Growth Factor 1 Antibody	1 vial
Chromogen Substrate	1 x 7ml
Insulin like Growth Factor 1 Microplate (12 x 8 well strips)	1 unit
Insulin like Growth Factor 1 Standard	1 vial
Sealing Tapes	1 x 3 units
Stop Solution	1 x 11ml

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.

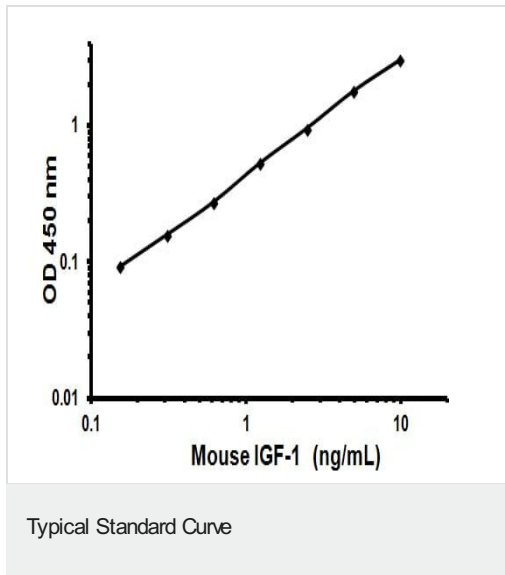
配列類似性

Belongs to the insulin family.

細胞内局在

Secreted.

画像



Representative Standard Curve using ab108874

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