

Rat IGF-2 ELISA Kit ab213903

2 References 1 图像

概述

产品名称 大鼠IGF-2 ELISA试剂盒

检测方法 Colorimetric

精确度 批次内

样品	n	Mean	SD	CV%
1	16	478pg/ml	20.55	4.3%
2	16	1536pg/ml	87.55	5.7%
3	16	2743pg/ml	126.2	4.6%

批次间

样品	n	Mean	SD	CV%
1	24	453pg/ml	30.8	6.8%
2	24	1737pg/ml	140.7	8.1%
3	24	2968pg/ml	222.6	7.5%

样品类型 Cell culture supernatant, Serum, Cell Lysate, Hep Plasma, EDTA Plasma, Tissue Homogenate

检测类型 Sandwich (quantitative)

灵敏度 < 5 pg/ml

范围 62.5 pg/ml - 4000 pg/ml

检测时间 3h 30m

实验步骤 Multiple steps standard assay

种属反应性 与反应: Rat

预测可用于: Hamster 

产品概述 The Rat IGF-2 Enzyme-Linked Immunosorbent Assay (ELISA) kit (ab213903) is designed for the quantitative measurement of Rat IGF-2 in cell culture supernatants, cell lysates, tissue homogenates, serum and plasma (heparin, EDTA).

The ELISA kit is based on standard sandwich enzyme-linked immune-sorbent assay technology. A monoclonal antibody from mouse specific for IGF-2 has been pre-coated onto 96-well plates. Standards and test samples are added to the wells, a biotinylated detection polyclonal antibody from goat specific for IGF-2 is added subsequently and then followed by washing with PBS or TBS buffer. Avidin-Biotin-Peroxidase Complex is added and unbound conjugates are washed away with PBS or TBS buffer. HRP substrate TMB is used to visualize HRP enzymatic reaction. TMB is catalyzed by HRP to produce a blue color product that changed into yellow after adding acidic stop solution. The density of yellow is proportional to the Rat IGF-2 amount of sample captured in plate.

## 说明

Insulin-like growth factor II (IGF-2) is also known as somatomedin A. IGF-2 is a member of the insulin family of polypeptide growth factors that is involved in development and growth. It is paternally expressed in the fetus and placenta. IGF-2 is a mitogen for many cell types and an important modulator of muscle growth and differentiation. IGF-2 gene is prevalently expressed during prenatal development and its gene activity is regulated by genomic imprinting, in that the allele inherited from the father is active and the allele inherited from the mother is inactive in most normal tissues. IGF-2 appears to be induced by placental lactogen during prenatal development. It is a mediator of prolactin-induced alveologenesis; prolactin, IGF-2, and cyclin D1, all of which are overexpressed in breast cancers, are components of a developmental pathway in the mammary gland.

## 平台

Pre-coated microplate (12 x 8 well strips)

## 性能

### 存放说明

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

组件	标识符	1 x 96 tests
ABC Diluent Buffer	Blue Cap	1 x 12ml
Adhesive Plate Seal		4 units
Antibody Diluent Buffer	Green Cap	1 x 12ml
Anti-rat IGF-2 coated Microplate (12 x 8 wells)		1 unit
Avidin-Biotin-Peroxidase Complex (ABC)		1 x 100µl
Biotinylated anti- Rat IGF-2 antibody		1 x 100µl
Lyophilized recombinant Rat IGF-2 standard		2 vials
Sample Diluent Buffer	Green Cap	1 x 30ml
TMB Color Developing Agent	Black Cap	1 x 10ml
TMB Stop Solution	Yellow Cap	1 x 10ml
Wash Buffer (25X)		1 x 20ml

## 功能

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.

#### 疾病相关

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.

#### 序列相似性

Belongs to the insulin family.

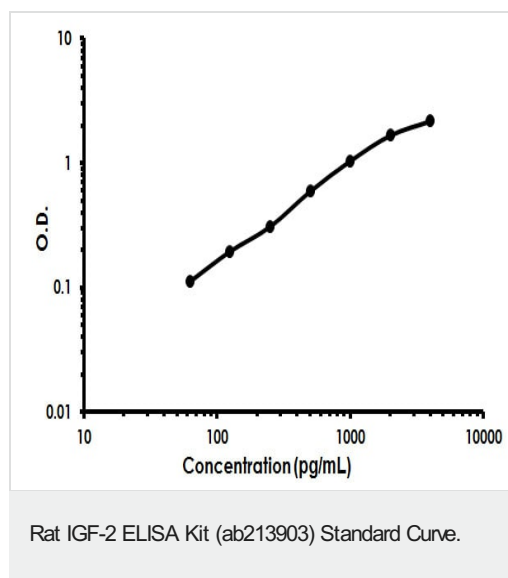
#### 翻译后修饰

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

#### 细胞定位

Secreted.

#### 图片



Rat IGF-2 ELISA Kit (ab213903) 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.cn/abpromise> or contact our technical team.

#### Terms and conditions

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