

Human Insulin like Growth Factor 1 ELISA Kit (IGF1) ab108873

★★★★★ 1 Abreviews 5 References 1 图像

概述

产品名称

检测方法

精确度

样品类型

检测类型

灵敏度

范围

回收率

检测时间

实验步骤

种属反应性

产品概述

人Insulin like Growth Factor 1 ELISA试剂盒(IGF1)

Colorimetric

批次内

样品	n	Mean	SD	CV%
Overall				6%

批次间

样品	n	Mean	SD	CV%
Overall				10.6%

Cell culture supernatant, Serum, Plasma

Sandwich (quantitative)

> 1.2 ng/ml

0.4 ng/ml - 24 ng/ml

= 98 %

5h 00m

Multiple steps standard assay

与反应: Human

Abcam's Insulin like Growth Factor 1 (IGF1) Human *in vitro* ELISA (Enzyme-Linked

Get results in 90 minutes with Human IGF1 ELISA Kit ([ab211651](#)) from our SimpleStep ELISA® range.

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

平台 Microplate

性能

存放说明 Store at -20°C. Please refer to protocols.

组件	1 x 96 tests
100X Streptavidin-Peroxidase Conjugate	1 x 80µl
10X Diluent N Concentrate	1 x 30ml
1X Plasma/Serum Pretreatment Buffer	1 x 7.5ml
1X Standard Diluent	1 x 2ml
20X Wash Buffer Concentrate	2 x 30ml
50X Biotinylated Human Insulin like Growth Factor 1 Antibody	1 x 120µl
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	3 units
Stop Solution	1 x 11ml

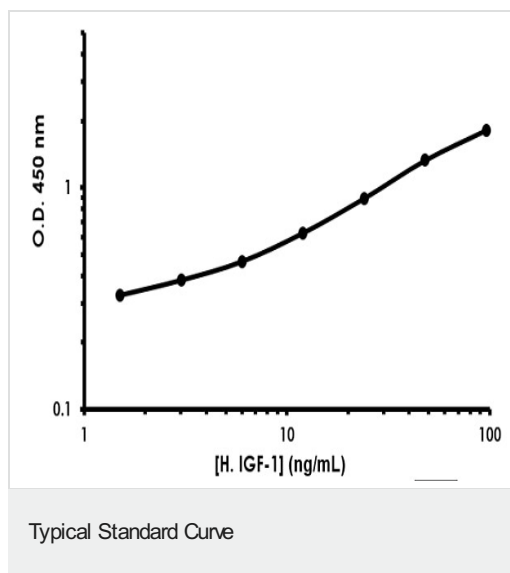
功能 The insulin-like growth factors, isolated from plasma, are structurally and functionally related to insulin but have a much higher growth-promoting activity. May be a physiological regulator of [1-14C]-2-deoxy-D-glucose (2DG) transport and glycogen synthesis in osteoblasts. Stimulates glucose transport 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.

疾病相关 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.

序列相似性 Belongs to the insulin family.

细胞定位 Secreted.

图片



Representative Standard Curve using ab108873.

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