

Human ARP4 ELISA Kit (Angiopoietin-Like 4) ab99974

9 References 2 图像

概述

产品名称	人ARP4 ELISA试剂盒(Angiopoietin-Like 4)
检测方法	Colorimetric
样品类型	Cell culture supernatant, Serum, Plasma
检测类型	Sandwich (quantitative)
灵敏度	< 20 pg/ml
范围	27.4 pg/ml - 20000 pg/ml
回收率	100.4 %

特定样本回收率

样品类型	平均%	范围
Cell culture supernatant	117.8	67% - 137%
Serum	95.38	85% - 104%
Plasma	88.04	74% - 103%

实验步骤	Multiple steps standard assay
种属反应性	与反应: Human
产品概述	<p>Abcam's ARP4 (Angiopoietin-like 4) Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an <i>in vitro</i> enzyme-linked immunosorbent assay for the quantitative measurement of Human ARP4 in serum, plasma, and cell culture supernatants.</p> <p>This assay employs an antibody specific for Human ARP4 coated on a 96-well plate. Standards and samples are pipetted into the wells and ARP4 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human ARP4 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 ARP4 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.</p>
说明	Optimisation may be required with urine samples.
平台	Microplate

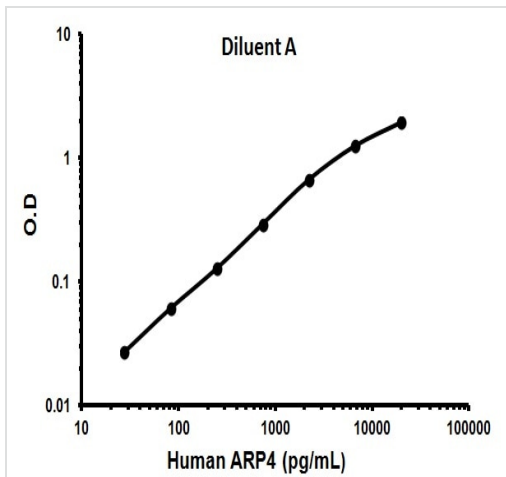
性能

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

组件	1 x 96 tests
20X Wash Buffer	1 x 25ml
280X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent B	1 x 15ml
ARP4 Microplate (12 x 8 well strips)	1 unit
Assay Diluent A	1 x 30ml
Biotinylated anti-Human ARP4	2 vials
Recombinant ARP4 Standard Human (Lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

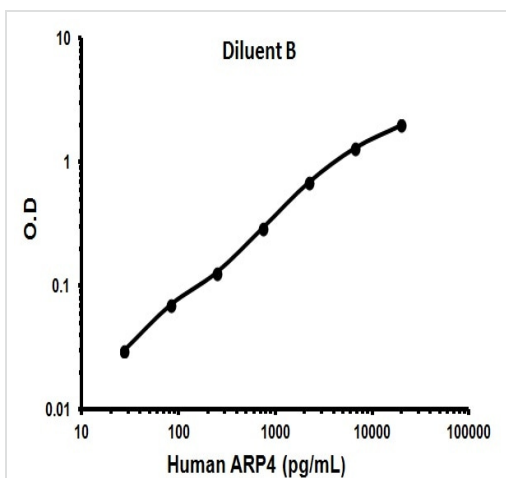
相关性 ARP4 is a fasting-induced adipose factor. It was identified as an adipocytokine up-regulated by fasting, by peroxisome proliferator-activated receptor agonists, and by hypoxia. At the protein level, in human and mouse blood plasma, ARP4 was found to be present both as a native protein and in a truncated form. Differentiation of mouse 3T3-L1 adipocytes was associated with the production of truncated ARP4, whereas in human white adipose tissue and SGBS adipocytes, only the native ARP4 could be detected. Interestingly, the truncated ARP4 was produced by human liver. Experimental data suggest that ARP4 is mainly presented in human blood plasma in a truncated form (ARP4-S2), whose level is increased by fenofibrate treatment. Levels of both truncated and native ARP4 showed marked inter individual variation but were not associated with body mass index and were not influenced by prolonged semistarvation.

图片



Representative standard curve using ab99974.

Typical Standard Curve



Representative standard curve using ab99974.

Typical 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