

Mouse Epiregulin ELISA Kit (EREG) ab100682

2 图像

概述

产品名称	小鼠Epiregulin ELISA试剂盒(EREG)
检测方法	Colorimetric
样品类型	Cell culture supernatant, Plasma
检测类型	Sandwich (quantitative)
灵敏度	< 10 pg/ml
范围	8.23 pg/ml - 6000 pg/ml
回收率	83 %

特定样本回收率

样品类型	平均%	范围
Cell culture supernatant	102.6	94% - 110%
Plasma	73.94	67% - 90%

实验步骤	Multiple steps standard assay
种属反应性	与反应: Mouse 不与反应: Human

**产品概述**

Abcam's Epiregulin Mouse ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Mouse Epiregulin in plasma and cell culture supernatants.

This assay employs an antibody specific for Mouse Epiregulin coated on a 96- well plate. Standards and samples are pipetted into the wells and Epiregulin present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Mouse Epiregulin 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 Epiregulin bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

We have not been able to detect endogenous Mouse Epiregulin in normal serum with ab100682, only in serum spiked with Mouse Epiregulin.

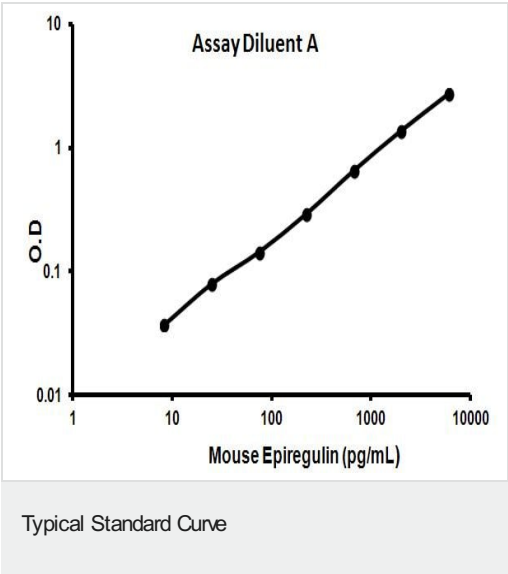
性能

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

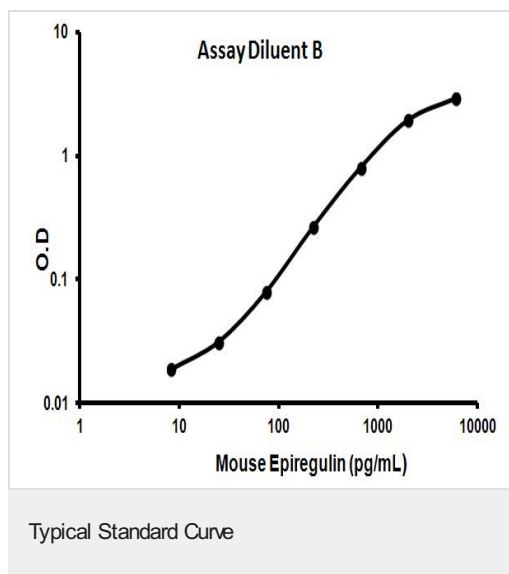
组件	1 x 96 tests
20X Wash Buffer	1 x 25ml
300X HRP-Streptavidin Concentrate	1 x 200μl
5X Assay Diluent B	1 x 15ml
Assay Diluent A	1 x 30ml
Biotinylated anti-Mouse Epiregulin	2 vials
Epiregulin Microplate (12 x 8 wells)	1 unit
Recombinant Mouse Epiregulin Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

相关性Epiregulin is a member of the EGF-related growth factor family. It is a mitogen and has numerous roles including as an insulinotropic factor in pancreatic beta cell lines. It may also be involved in metastasis.

图片



Representative Standard Curve using ab100682



Representative Standard Curve using ab100682

**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