

Mouse Epiregulin ELISA Kit (EREG) ab100682

2 Images

Overview

Product name	Mouse Epiregulin ELISA Kit (EREG)
Detection method	Colorimetric
Sample type	Cell culture supernatant, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	< 10 pg/ml
Range	8.23 pg/ml - 6000 pg/ml
Recovery	83 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	102.6	94% - 110%
Plasma	73.94	67% - 90%

Assay duration	Multiple steps standard assay
Species reactivity	Reacts with: Mouse Does not react with: Human

Product overview

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.

Properties

Storage instructions

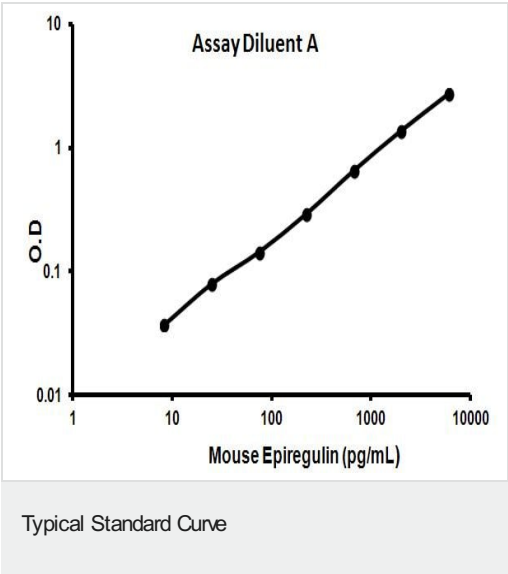
Store at -20°C. Please refer to protocols.

Components	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

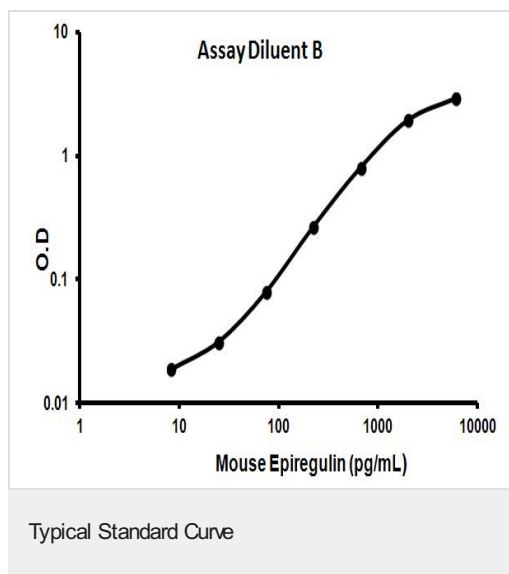
Relevance

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.

Images



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.com/abpromise> or contact our technical team.

Terms and conditions

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