

Human Amphiregulin ELISA Kit (AREG) ab99975

6 References [画像数 2](#)

製品の概要

製品名	Human Amphiregulin ELISA Kit (AREG)
検出方法	Colorimetric
サンプルの種類	Cell culture supernatant, Plasma
アッセイタイプ	Sandwich (quantitative)
検出感度	< 10 pg/ml
検出範囲	16.46 pg/ml - 4000 pg/ml
添加回収試験	83 %

特定サンプルでの回収試験

サンプルの種類	平均 %	測定範囲
Cell culture supernatant	100.4	89% - 111%
Plasma	72.7	64% - 87%

ステップ Multiple steps standard assay

種交差性 交差種: Human

製品の概要 Abcam's Amphiregulin (AREG) Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Human Amphiregulin in plasma, and cell culture supernatants. (Human Amphiregulin concentration is pretty low in normal plasma, it may not be detected in this assay).

This assay employs an antibody specific for Human Amphiregulin coated on a 96-well plate. Standards and samples are pipetted into the wells and Amphiregulin present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human Amphiregulin 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 Amphiregulin 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 Human Amphiregulin in normal serum with ab99975, only in serum spiked with Human Amphiregulin.

特記事項 Optimisation may be required with urine samples.

試験プラットフォーム Microplate

製品の特性

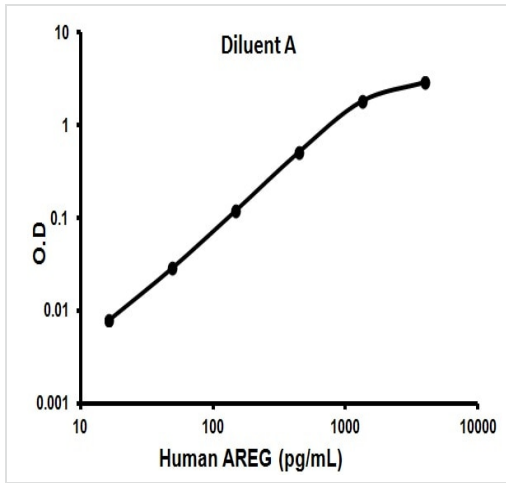
保存方法 Store at -20°C. Please refer to protocols.

内容	1 x 96 tests
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
500X HRP-Streptavidin Concentrate	1 x 200µl
Amphiregulin Microplate (12 x 8 well strips)	1 unit
Assay Diluent A	1 x 30ml
Biotinylated anti-Human Amphiregulin	2 vials
Recombinant Human Amphiregulin Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

関連性 The mature secreted form of amphiregulin is an 84-amino acid residue glycosylated polypeptide which is generated by proteolytic processing of a 252-amino acid transmembrane precursor. Seven different polypeptide ligands, which derive from distinct genes, are capable of binding to the extracellular domain of EGFR; these are EGF, TGF α , AR, HB-EGF, cripto-1, epiregulin, and betacellulin. All of these growth factors contain a characteristic EGF-like domain which is defined by 6 evenly spaced cysteine residues that generate 3 loops through the formation of disulfide bonds. Amphiregulin expression is induced by phorbol ester, estrogen, androgen, and other EGFR ligands. In vitro, amphiregulin functions as an autocrine growth stimulator to drive proliferation of colon carcinoma cells, normal and oncogene-transformed mammary epithelial cells, cervical carcinoma cells, prostate cancer cells, and keratinocytes. Amphiregulin is important in the development of human placenta and murine mammary gland. It is reportedly overexpressed in human cancers of breast, colon, stomach and pancreas.

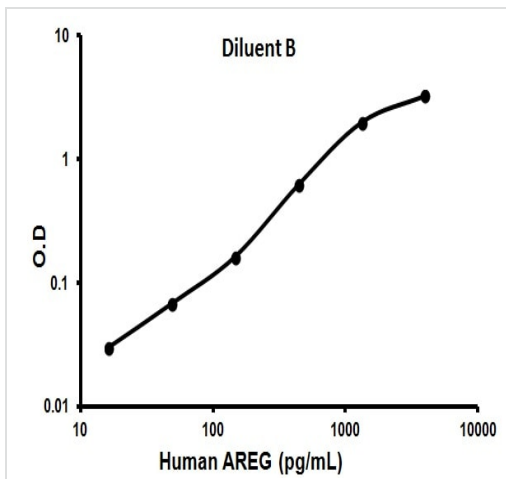
細胞内局在 Cell Membrane

画像



Representative Standard Curve using ab99975.

Typical Standard Curve



Representative Standard Curve using ab99975.

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.co.jp/abpromise> or contact our technical team.

Terms and conditions

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