

Mouse Amphiregulin ELISA Kit α b100668

3 References 1 Image

Overview

Product name	Mouse Amphiregulin ELISA Kit
Detection method	Colorimetric
Sample type	Cell culture supernatant, Serum, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	< 2 pg/ml
Range	1.1 pg/ml - 800 pg/ml
Recovery	100 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	85.25	76% - 95%
Serum	130.1	120% - 138%
Plasma	123.4	112% - 130%

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Mouse

Product overview Abcam’s Amphiregulin Mouse ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of mouse Amphiregulin in serum, plasma and cell culture supernatants.

This assay employs an antibody specific for mouse 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-mouse 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.

Platform Microplate

Properties

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

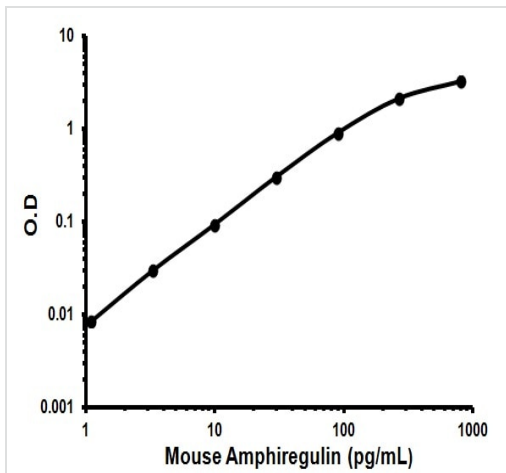
Components	1 x 96 tests
200X HRP-Streptavidin Concentrate	1 x 200µl
20X Wash Buffer	1 x 25ml
5X Assay Diluent	1 x 15ml
Amphiregulin Microplate (12 x 8 wells)	1 unit
Biotinylated anti-Mouse Amphiregulin	2 vials
Recombinant Mouse Amphiregulin Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step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

Relevance

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.

Cellular localization Cell Membrane

Images



Representative Standard Curve using ab100668.

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.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