

Product datasheet

Human Amphiregulin ELISA Kit ab222504

Recombinant SimpleStep ELISA

[5 Images](#)

Overview

Product name Human Amphiregulin ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

| Sample | n | Mean | SD | CV% |
|--------------|---|------|----|------|
| Amphiregulin | 3 | | | 1.8% |

Inter-assay

| Sample | n | Mean | SD | CV% |
|--------------|---|------|----|------|
| Amphiregulin | 5 | | | 4.1% |

Sample type

Cell culture supernatant, Serum, EDTA Plasma, Cit plasma

Assay type

Sandwich (quantitative)

Sensitivity

2.4 pg/ml

Range

18 pg/ml - 600 pg/ml

Recovery

Sample specific recovery

| Sample type | Average % | Range |
|--------------------|-----------|-------------|
| Serum | 106 | 84% - 117% |
| Cell culture media | 108 | 106% - 110% |
| EDTA Plasma | 98 | 91% - 104% |
| Cit plasma | 103 | 93% - 111% |

Assay time

1h 30m

Assay duration

One step assay

Species reactivity**Reacts with:** Human**Product overview**

Human Amphiregulin ELISA Kit (ab222504) is a single-wash 90 min sandwich ELISA designed for the quantitative measurement of Amphiregulin protein in cell culture supernatant, cit plasma, edta plasma, and serum. It uses our proprietary SimpleStep ELISA® technology. Quantitate Human Amphiregulin with 2.4 pg/ml sensitivity.

SimpleStep ELISA® technology employs capture antibodies conjugated to an affinity tag that is recognized by the monoclonal antibody used to coat our SimpleStep ELISA® plates. This approach to sandwich ELISA allows the formation of the antibody-analyte sandwich complex in a single step, significantly reducing assay time. See the SimpleStep ELISA® protocol summary in the image section for further details. Our SimpleStep ELISA® technology provides several benefits:

- Single-wash protocol reduces assay time to 90 minutes or less
- High sensitivity, specificity and reproducibility from superior antibodies
- Fully validated in biological samples
- 96-wells plate breakable into 12 x 8 wells strips

A 384-well SimpleStep ELISA® microplate ([ab203359](#)) is available to use as an alternative to the 96-well microplate provided with SimpleStep ELISA® kits.

Notes

Amphiregulin is a member of the epidermal growth factor family encoded by the AREG gene. Amphiregulin is expressed as a 252- amino acid transmembrane glycoprotein, which contains an 81-amino acid N-terminal pro-peptide, a heparin binding domain, a EGF like domain, a 23- amino acid transmembrane domain, and a 31- amino acid cytoplasmic domain. The amphiregulin assay utilizes antibodies raised against the 101-187 amino acid region of the amphiregulin protein. Amphiregulin is an autocrine growth factor, a mitogen for several cell types, and is essential for mammary ductal development.

Platform

Pre-coated microplate (12 x 8 well strips)

Properties**Storage instructions**

Store at +4°C. Please refer to protocols.

| Components | 1 x 96 tests |
|---|--------------|
| 10X Human Amphiregulin Capture Antibody | 1 x 600µl |
| 10X Human Amphiregulin Detector Antibody | 1 x 600µl |
| 10X Wash Buffer PT (ab206977) | 1 x 20ml |
| Antibody Diluent CPI - HAMA Blocker (ab193969) | 1 x 6ml |
| Human Amphiregulin Lyophilized Recombinant Protein | 2 vials |
| Plate Seals | 1 unit |
| SimpleStep Pre-Coated 96-Well Microplate (ab206978) | 1 unit |

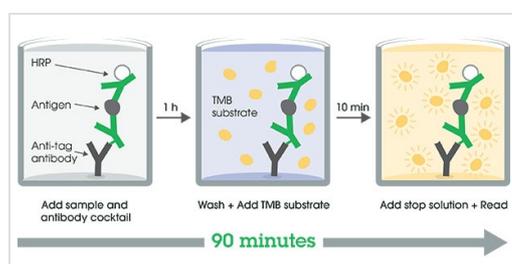
| Components | 1 x 96 tests |
|------------------------------|--------------|
| Sample Diluent NS (ab193972) | 1 x 50ml |
| Stop Solution | 1 x 12ml |
| TMB Development Solution | 1 x 12ml |

Function Bifunctional growth-modulating glycoprotein. Inhibits growth of several human carcinoma cells in culture and stimulates proliferation of human fibroblasts and certain other tumor cells.

Sequence similarities Belongs to the amphiregulin family.
Contains 1 EGF-like domain.

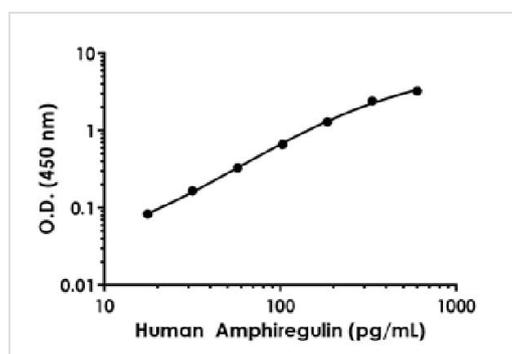
Cellular localization Membrane.

Images



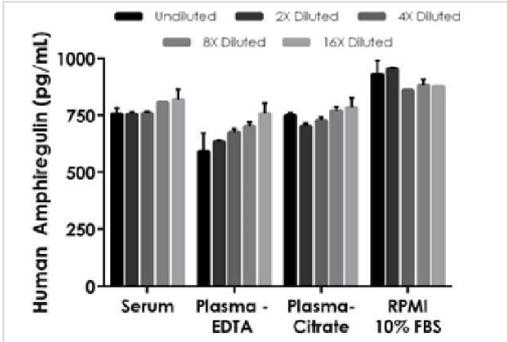
Other - Human Amphiregulin ELISA Kit (ab222504)

SimpleStep ELISA technology allows the formation of the antibody-antigen complex in one single step, reducing assay time to 90 minutes. Add samples or standards and antibody mix to wells all at once, incubate, wash, and add your final substrate. See protocol for a detailed step-by-step guide.



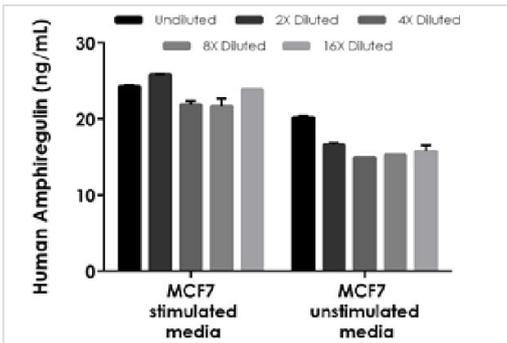
Example of human Amphiregulin standard curve in Sample Diluent NS.

The Amphiregulin standard curve was prepared as described in Section 10. Raw data values are shown in the table. Background-subtracted data values (mean +/- SD) are graphed.



Interpolated concentrations of spiked Amphiregulin in human serum, plasma and cell culture supernatant samples.

The concentrations of Amphiregulin were measured in duplicates, interpolated from the Amphiregulin standard curves and corrected for sample dilution. Undiluted samples are as follows: serum 50%, plasma (EDTA) 50%, plasma (citrate) 50%, RPMI media 50%. The interpolated dilution factor corrected values are plotted (mean \pm SD, n=2).



Interpolated concentrations of native Amphiregulin in human MCF7 cell culture supernatant.

Human MCF7s were stimulated with 2% PHA for 3 days before the supernatant was collected and assayed. The concentrations of Amphiregulin were measured in duplicate and interpolated from the Amphiregulin standard curve and corrected for sample dilution. The interpolated dilution factor corrected values are plotted (mean \pm SD, n=2). The mean Amphiregulin concentration was determined to be 23.7 ng/mL in stimulated and 16 ng/mL in unstimulated MCF7 cell culture supernatant.

Powered by
recombinant antibodies



Research with confidence
Consistent and reproducible results



Long-term and scalable supply
Recombinant technology



Success from the first experiment
Confirmed specificity



Ethical standards compliant
Animal-free production

Sandwich ELISA - Human Amphiregulin ELISA Kit
(ab222504)

To learn more about the advantages of recombinant antibodies see [here](#).

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