abcam

Product datasheet

Human Epiregulin ELISA Kit ab213775

1 Image

Overview

Product name Human Epiregulin ELISA Kit

Detection methodColorimetric

Precision Intra-assay

Sample	n	Mean	SD	CV%
1	16	13pg/ml	0.78	= 6%
2	16	57pg/ml	4.21	= 7.4%
3	16	113pg/ml	6.1	= 4.6%

Inter-assay

Sample	n	Mean	SD	CV%
1	24	12pg/ml	0.94	= 7.9%
2	24	56pg/ml	4.48	= 8%
3	24	105pg/ml	6.4	= 6.1%

Sample type Cell culture supernatant, Serum, Cell Lysate, Hep Plasma, EDTA Plasma

Assay type Sandwich (quantitative)

Sensitivity < 10 pg/ml

Range 3.9 pg/ml - 250 pg/ml

Assay time 3h 30m

Assay duration Multiple steps standard assay

Species reactivity Reacts with: Human

Product overview The Human Epiregulin Enzyme-Linked Immunosorbent Assay (ELISA) kit (ab213775) is designed

for the quantitative measurement of Human Epiregulin in cell culture supernatants, cell lysates,

serum and plasma (heparin, EDTA).

The ELISA kit is based on standard sandwich enzyme-linked immune-sorbent assay technology.

A monoclonal antibody from mouse specific for Epiregulin has been pre-coated onto 96-well plates. Standards (*E.coli*; V60-L108) and test samples are added to the wells, a biotinylated detection polyclonal antibody from goat specific for Epiregulin is added subsequently and then followed by washing with PBS or TBS buffer. Avidin-Biotin-Peroxidase Complex is added and unbound conjugates are washed away with PBS or TBS buffer. HRP substrate TMB is used to visualize HRP enzymatic reaction. TMB is catalyzed by HRP to produce a blue color product that changed into yellow after adding acidic stop solution. The density of yellow is proportional to the Human Epiregulin amount of sample captured in plate.

Notes

Epiregulin (EPR) is a protein that in humans is encoded by the EREG gene. It is a member of the epidermal growth factor family. Epiregulin can function as a ligand of epidermal growth factor receptor (EGFR), as well as a ligand of most members of the ERBB (v-erb-b2 oncogene homolog) family of tyrosine-kinase receptors. The secondary structure at the C-terminus epiregulin is different from other epidermal growth factor family ligands because of the lack of hydrogen bonds. The structural difference at the C-terminus may provide an explanation for the reduced binding affinity of epiregulin to the ERBB receptors.

Platform

Pre-coated microplate (12 x 8 well strips)

Properties

Storage instructions

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

Components	Identifier	1 x 96 tests
ABC Diluent Buffer	Blue Cap	1 x 12ml
Adhesive Plate Seal		4 units
Antibody Diluent Buffer	Green Cap	1 x 12ml
Anti-Human Epiregulin coated Microplate (12 x 8 wells)		1 unit
Avidin-Biotin-Peroxidase Complex (ABC)		1 x 100µl
Biotinylated anti- Human Epiregulin antibody		1 x 100µl
Lyophilized recombinant Human Epiregulin Standard		2 vials
Sample Diluent Buffer	Green Cap	1 x 30ml
TMB Color Developing Agent	Black Cap	1 x 10ml
TMB Stop Solution	Yellow Cap	1 x 10ml
Wash Buffer (25X)		1 x 20ml

Function May be a mediator of localized cell proliferation. As a mitogen it may stimulate cell proliferation

and/or angiogenesis.

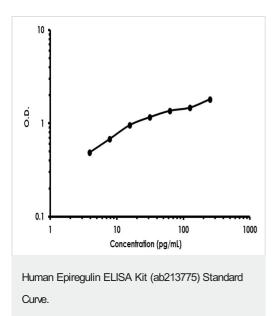
Tissue specificity In normal adults, expressed predominantly in the placenta and peripheral blood leukocytes. High

levels were detected in carcinomas of the bladder, lung, kidney and colon.

Sequence similarities Contains 1 EGF-like domain.

2

Images



Human Epiregulin ELISA Kit (ab213775) 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