

Human BAFF-R ELISA Kit ab213839

1 Image

Overview

Product name Human BAFF-R ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
1	16	147pg/ml	10.29	= 7%
2	16	683pg/ml	49.85	= 7.3%
3	16	2750pg/ml	206.25	= 7.5%

Inter-assay

Sample	n	Mean	SD	CV%
1	24	156pg/ml	13.88	= 8.9%
2	24	700pg/ml	62.3	= 8.9%
3	24	2531pg/ml	227.05	= 8.9%

Sample type

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

Assay type

Sandwich (quantitative)

Sensitivity

< 10 pg/ml

Range

78 pg/ml - 5000 pg/ml

Assay time

3h 30m

Assay duration

Multiple steps standard assay

Species reactivity

Reacts with: Human

Product overview

The Human BAFF-R Enzyme-Linked Immunosorbent Assay (ELISA) kit (ab213839) is designed for the quantitative measurement of Human TNFRSF13C 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 goat polyclonal antibody specific for BAFF-R has been pre-coated onto 96-well plates. Standards (Expression system for standard: NSO; Immunogen sequence: S7-A71) and test samples are added to the wells, a biotinylated detection polyclonal antibody from goat specific for BAFF-R 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 BAFF-R amount of sample captured in plate.

Notes BAFF-R (also known as Tumor necrosis factor receptor superfamily member 13C or B-cell-activating factor receptor) is a protein that in humans is encoded by the TNFRSF13C gene. By homology to a BAC clone, the BAFFR gene was mapped to chromosome 22q13.1-q13.31. B-cell activating factor (BAFF) enhances B-cell survival *in vitro* and is a regulator of the peripheral B-cell population. The protein encoded by this gene is a receptor for BAFF and is a type III transmembrane protein containing a single extracellular phenylalanine-rich domain. It is thought that this receptor is the principal receptor required for BAFF-mediated mature B-cell survival.

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 BAFF-R coated Microplate (12 x 8 wells)		1 unit
Avidin-Biotin-Peroxidase Complex (ABC)		1 x 100µl
Biotinylated anti-Human BAFF-R antibody		1 x 100µl
Lyophilized Recombinant Human BAFF-R 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 B-cell receptor specific for TNFSF13B/TALL1/BAFF/BLyS. Promotes the survival of mature B-cells and the B-cell response.

Tissue specificity Highly expressed in spleen and lymph node, and in resting B-cells. Detected at lower levels in activated B-cells, resting CD4+ T-cells, in thymus and peripheral blood leukocytes.

Involvement in disease Defects in TNFRSF13C are the cause of immunodeficiency common variable type 4 (CVID4)

[MIM:613494]; also called antibody deficiency due to BAFFR defect. CVID4 is a primary immunodeficiency characterized by antibody deficiency, hypogammaglobulinemia, recurrent bacterial infections and an inability to mount an antibody response to antigen. The defect results from a failure of B-cell differentiation and impaired secretion of immunoglobulins; the numbers of circulating B cells is usually in the normal range, but can be low.

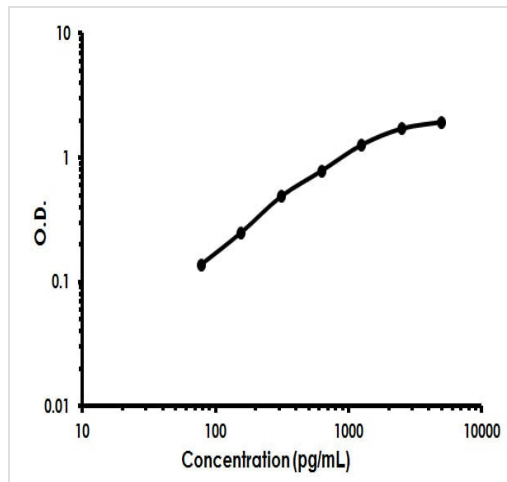
Sequence similarities

Contains 1 TNFR-Cys repeat.

Cellular localization

Membrane.

Images



Human TNFRSF13C ELISA Kit (BAFFR) (ab213839) Standard Curve.

Human TNFRSF13C ELISA Kit (BAFFR) (ab213839) 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