



Product datasheet

Human BAFF ELISA Kit (TNFSF13B) ab188391

SimpleStep ELISA

3 References 5 图像

概述

产品名称 人BAFF ELISA试剂盒(TNFSF13B)

检测方法 Colorimetric

精确度					批次内
	样品	n	Mean	SD	CV%
	Overall	8			8.3%

					批次间
	样品	n	Mean	SD	CV%
	Overall	3			8.8%

样品类型 Serum, Hep Plasma, EDTA Plasma, Cit plasma

检测类型 Sandwich (quantitative)

灵敏度 12.7 pg/ml

范围 0.08 ng/ml - 5 ng/ml

回收率			特定样本回收率
	样品类型	平均%	范围
	Serum	111	105% - 115%
	Cell culture media	108	106% - 110%
	Hep Plasma	104	101% - 108%
	EDTA Plasma	105	98% - 115%
	Cit plasma	106	100% - 111%

检测时间 1h 30m

实验步骤 One step assay

产品概述

Human BAFF ELISA Kit (TNFSF13B) (ab188391) is a single-wash 90 min sandwich ELISA designed for the quantitative measurement of BAFF (TNFSF13B) protein in cit plasma, edta plasma, hep plasma, and serum. It uses our proprietary SimpleStep ELISA® technology. Quantitate Human BAFF (TNFSF13B) with 12.7 pg/ml sensitivity.

- 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

说明

Inadequate levels of BAFF will fail to activate B-cells to produce enough immunoglobulin and will lead to immunodeficiency. Excessive level of BAFF causes abnormally high antibody production, results in systemic lupus erythmatosis, rheumatoid arthritis, and many other autoimmune diseases.

平台

Microplate

性能

存放说明

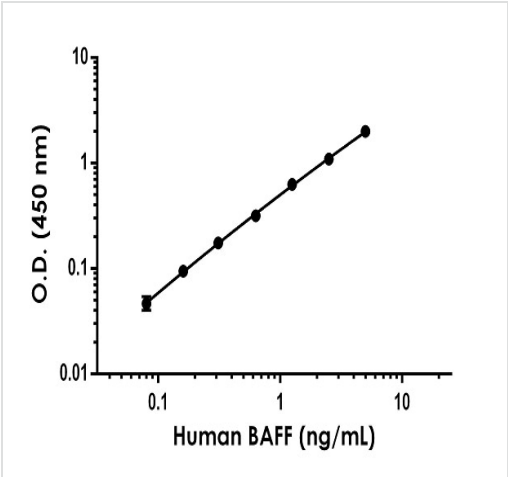
Store at +4°C. Please refer to protocols.

组件	1 x 96 tests
10X Human BAFF Capture Antibody	1 x 600μl
10X Human BAFF Detector Antibody	1 x 600μl
10X Wash Buffer PT (ab206977)	1 x 20ml
Antibody Diluent 4BI	1 x 6ml

组件	1 x 96 tests
Human BAFF Lyophilized Recombinant Protein	2 vials
Plate Seals	1 unit
Sample Diluent 25BS	1 x 20ml
Sample Diluent NS (ab193972)	1 x 12ml
SimpleStep Pre-Coated 96-Well Microplate (ab206978)	1 unit
Stop Solution	1 x 12ml
TMB Development Solution	1 x 12ml

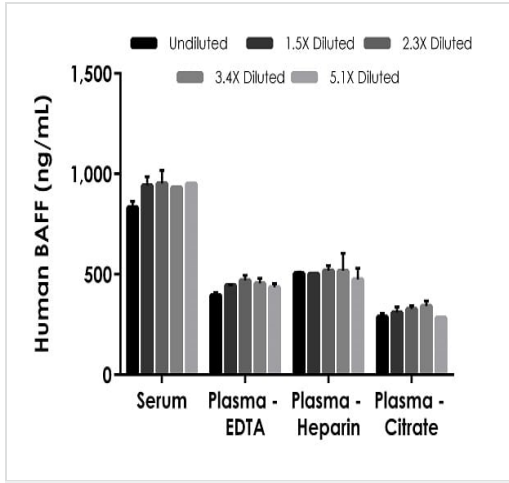
功能	Cytokine that binds to TNFRSF13B/TACI and TNFRSF17/BCMA. TNFSF13/APRIL binds to the same 2 receptors. Together, they form a 2 ligands -2 receptors pathway involved in the stimulation of B-and T-cell function and the regulation of humoral immunity. A third B-cell specific BAFF-receptor (BAFFR/BR3) promotes the survival of mature B-cells and the B-cell response.
组织特异性	Abundantly expressed in peripheral blood Leukocytes and is specifically expressed in monocytes and macrophages. Also found in the spleen, lymph node, bone marrow, T-cells and dendritic cells. A lower expression seen in placenta, heart, lung, fetal liver, thymus, and pancreas.
序列相似性	Belongs to the tumor necrosis factor family.
翻译后修饰	The soluble form derives from the membrane form by proteolytic processing. N-glycosylated.
细胞定位	Secreted and Cell membrane.

图片



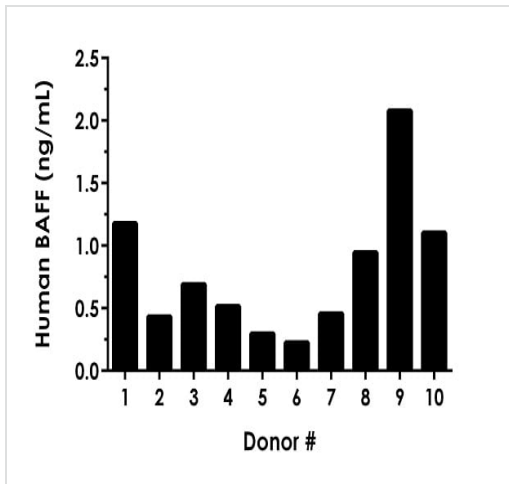
Example of BAFF standard curve.

Background-subtracted data values (mean +/- SD) are graphed.



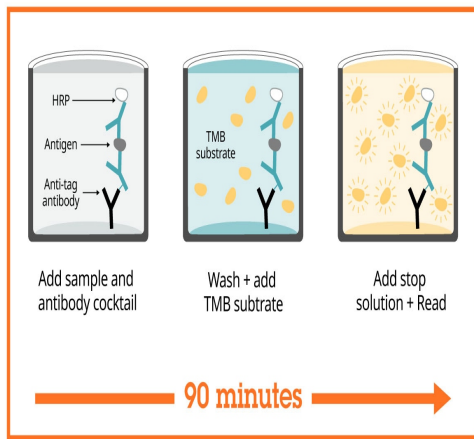
Background-subtracted data values (mean +/- SD, n = 2) are graphed.

Titration of Human Plasma Samples within the working range of the assay.



Quantification of BAFF interpolated from standard curve and multiplied by dilution factor.

Titration of individual donor serum samples to within working range of the assay.



Sandwich ELISA - Human BAFF ELISA Kit
(TNFSF13B) (ab188391)

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.

Get more done with SimpleStep ELISA



Easy to use
Single-wash 90-minute protocol



Flexible
Matched antibody pairs available



Precision antibodies
High sensitivity, specificity and reproducibility



Scalable
Now in 10-pack and 384-well formats

Sandwich ELISA - Human BAFF ELISA Kit
(TNFSF13B) (ab188391)

To learn more about the advantages of SimpleStep ELISA® kits 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.cn/abpromise> or contact our technical team.

Terms and conditions

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