

Human BAFF ELISA Kit (TNFSF13B) ab119579

2 References 1 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%
	1	16	127pg/ml	5.46	= 4.3%
	2	16	522pg/ml	38.62	= 7.4%
	3	16	1229pg/ml	79.88	= 6.5%
	批次间				
	样品	n	Mean	SD	CV%
	1	24	124pg/ml	5.7	= 4.6%
	2	24	548pg/ml	47.67	= 8.7%
	3	24	1268pg/ml	103.97	= 8.2%
样品类型	Cell culture supernatant, Serum, Plasma, Cell Lysate				
检测类型	Sandwich (quantitative)				
灵敏度	< 2 pg/ml				
范围	62.5 pg/ml - 4000 pg/ml				
实验步骤	Multiple steps standard assay				
种属反应性	与反应: Human				
产品概述	Abcam's Human BAFF (TNFSF13B) <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate quantitative measurement of Human BAFF (TNFSF13B) in cell culture supernatants, serum and plasma (heparin, EDTA).				

A BAFF (TNFSF13B) specific mouse monoclonal antibody has been precoated onto 96-well plates. Standards and test samples are added to the wells and incubated. A biotinylated

detection polyclonal antibody from goat, specific for BAFF (TNFSF13B) is then added 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. TMB is then used to visualize the HRP enzymatic reaction. TMB is catalyzed by HRP to produce a blue color product that changes into yellow after adding acidic stop solution. The density of yellow coloration is directly proportional to the Human BAFF (TNFSF13B) amount of sample captured in plate.

Get higher sensitivity in only 90 minutes with Human BAFF ELISA Kit (TNFSF13B) (**ab188391**) from our SimpleStep ELISA® range.

**平台** Microplate

## 性能

**存放说明** Store at -20°C. Please refer to protocols.

组件	标识符	1 x 96 tests	1 x 96 tests
ABC Diluent Buffer	Blue Cap	1 x 12ml	1 x 12ml
Antibody Diluent Buffer	Green Cap	1 x 12ml	1 x 12ml
Anti-Human BAFF Antibody Microplate (12 x 8 wells)		1 unit	1 unit
Avidin-Biotin-Peroxidase Complex (ABC)		1 x 100µl	1 x 100µl
Biotinylated anti-Human BAFF antibody		1 x 100µl	1 x 100µl
Lyophilized recombinant Human BAFF standard		2 x 10ng	2 x 10ng
Plate Seal		4 units	4 units
Sample Diluent Buffer	Green Cap	1 x 30ml	1 x 30ml
TMB Color Developing Agent	Amber Bottle	1 x 10ml	1 x 10ml
TMB Stop Solution	Yellow Cap	1 x 10ml	1 x 10ml
Wash Buffer (25X)		1 x 20ml	1 x 20ml

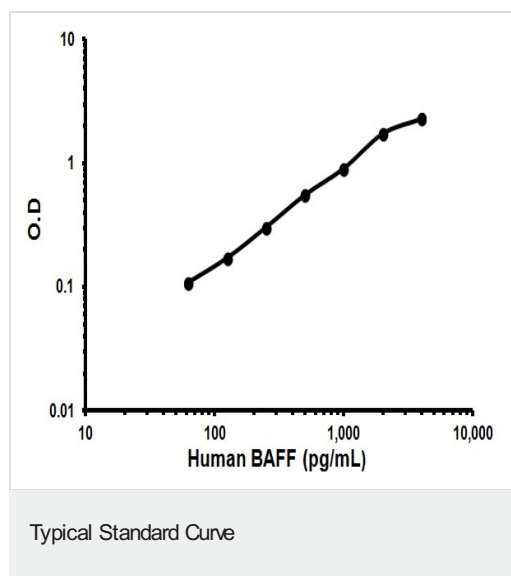
**功能** Cytokine that binds to TNFRSF13B/TACI and TNFRSF17/BCMA. TNFSF13/APRIL binds to the same 2 receptors. Together, they form a 2 ligands -2 receptors pathway involved in the stimulation of B-and T-cell function and the regulation of humoral immunity. A third B-cell specific BAFF-receptor (BAFFR/BR3) 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.



Representative Standard Curve using ab119579.

**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