

Human BAFF ELISA Kit (TNFSF13B) ab119579

2 References 画像数 1

製品の概要

製品名	Human BAFF ELISA Kit (TNFSF13B)			
検出方法	Colorimetric			
再現性	Intra-Assay (同時再現性)			
	サンプル	N	平均値	SD
	1	16	127pg/ml	5.46
	2	16	522pg/ml	38.62
	3	16	1229pg/ml	79.88
				CV%
				= 4.3%
				= 7.4%
				= 6.5%
	Inter-Assay (日差再現性)			
	サンプル	N	平均値	SD
	1	24	124pg/ml	5.7
	2	24	548pg/ml	47.67
	3	24	1268pg/ml	103.97
				CV%
				= 4.6%
				= 8.7%
				= 8.2%
サンプルの種類	Cell culture supernatant, Serum, Plasma, Cell Lysate			
アッセイタイプ	Sandwich (quantitative)			
検出感度	< 2 pg/ml			
検出範囲	62.5 pg/ml - 4000 pg/ml			
ステップ	Multiple steps standard assay			
種交差性	交差種: Human			
製品の概要	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.

#### 試験プラットフォーム

Microplate

#### 製品の特長

#### 保存方法

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

内容	ラベル	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

#### 機能

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.

#### 組織特異性

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.

#### 配列類似性

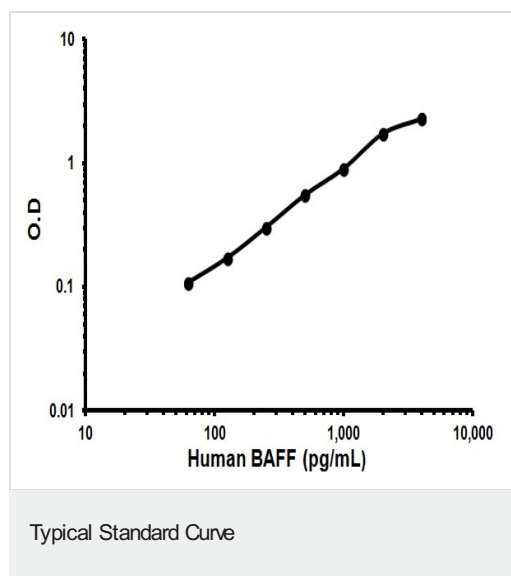
Belongs to the tumor necrosis factor family.

#### 翻訳後修飾

The soluble form derives from the membrane form by proteolytic processing. N-glycosylated.

#### 細胞内局在

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.co.jp/abpromise> or contact our technical team.

### Terms and conditions

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