

# BAFF soluble (mouse), detection set

BAFF (B cell activation factor of the TNF family, also known as BLyS or TALL1) is a cytokine expressed predominantly by cells of the immune system such as neutrophils, monocytes, macrophages, dendritic cells, follicular dendritic cells, activated T cells and some malignant B cells. BAFF binds three distinct receptors (BAFF-R, TACI and BCMA) expressed predominantly on B cells, although activated T cells also express BAFF-R. BAFF is a master regulator of peripheral B cell survival, and also acts in processes such as immunoglobulin isotype switch and B cell co-stimulation. Besides its major role in B cell biology, BAFF co-stimulates activated T cells. Deregulated expression of this membrane-bound protein, which can readily be released in a soluble form, leads to autoimmune disorders in mice. In humans, elevated levels of soluble BAFF have been detected in the serum of patients with various autoimmune diseases.

Citations: 5

[View Online »](#)

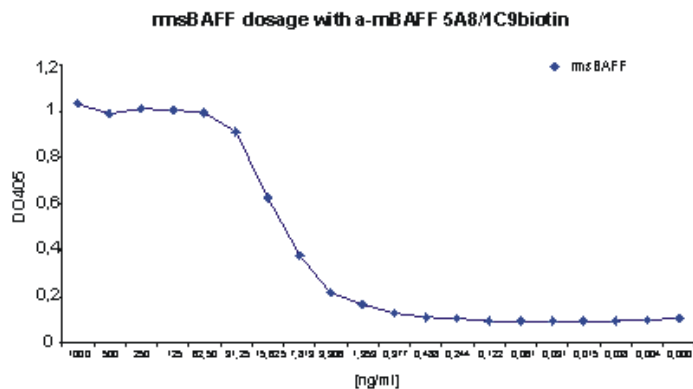
## Ordering Information

[Order Online »](#)

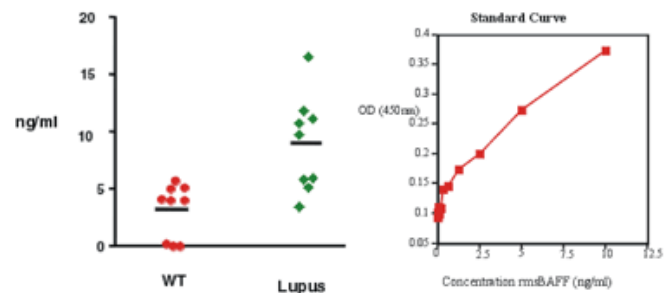
APO-54N-013-KI01	4x96 wells
------------------	------------

## Manuals, SDS & CofA

[View Online »](#)



**Figure:** Standard curve: coated with MAb to BAFF (mouse) (5A8) (Prod. No. [ALX-804-158](#)) and detected with MAb to BAFF (mouse) (1C9) (Biotin) (Prod. No. ALX-804-157B) followed by SAP-HRP. **Standard Curve Method:** A 96-well ELISA plate was coated O/N at RT with 100µl/well of a 5µg/ml solution of MAb to BAFF (mouse) (5A8) (Prod. No. ALX-804-158) diluted in PBS



**Figure:** left) Serum levels of BAFF (mouse) in wild-type (red) and lupus-prone mice (green) measured with the matched pair MAbs 5A8 and 1C9. right) Standard curve: coated with 5A8 (5µg/ml) and detected with 1C9 (Biotin) followed by SAP-HRP.

**Standard Curve Method:** A 96-well ELISA plate was coated O/N at 4°C with 100µl/well of a 5µg/ml solution of MAb to BAFF (mouse) (5A8) (Prod. No. ALX-804-158 ) diluted in 50mM sodium bicarbonate buffer (pH 9.6). After two washes

## Handling & Storage

<b>Use/Stability</b>	As indicated on product label or CoA when stored as recommended.
<b>Handling</b>	For long term storage keep the protein standard at -20°C. Avoid freeze/thaw cycles.
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Blue Ice

## Regulatory Status

RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	BLyS, TALL-1, THANK, zTNF4, TNFSF 13B/20, CD257
<b>Application</b>	Colorimetric detection, ELISA
<b>Application Notes</b>	For the quantitative determination of soluble mouse BAFF from biological fluids (serum and cell culture supernatant).
<b>Contents</b>	1 vial Standard (lyophilized) (STD) 1 vial Coating Antibody (COAT) 1 vial Detection Antibody (DET)
<b>Sensitivity</b>	0.3ng/ml (range 0 to 20ng/ml)
<b>UniProt ID</b>	Q9WU72
<b>Wavelength</b>	450 nm



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)