

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 »](#)

Manuals, SDS & CofA

[View Online »](#)

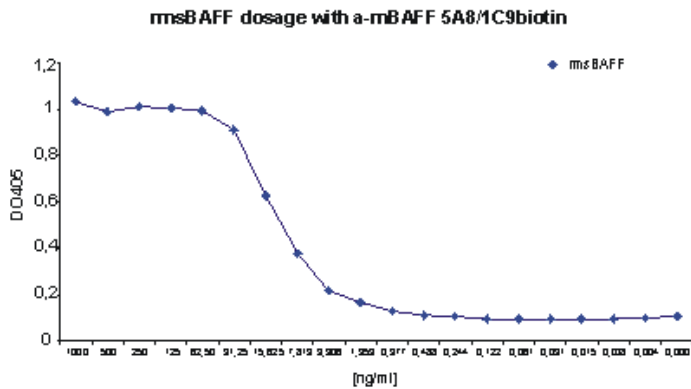


Figure: Standard curve: coated with MAb to BAFF (mouse) (5A8) (Prod. No. [ALX-804-158](#)) and detected with MAb to BAFF (mouse) (1C9) (Biotin) (Prod. No. ALX-804-157B) followed by SAP-HRP. **Standard Curve Method:** A 96-well ELISA plate was coated O/N at RT with 100µl/well of a 5µg/ml solution of MAb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

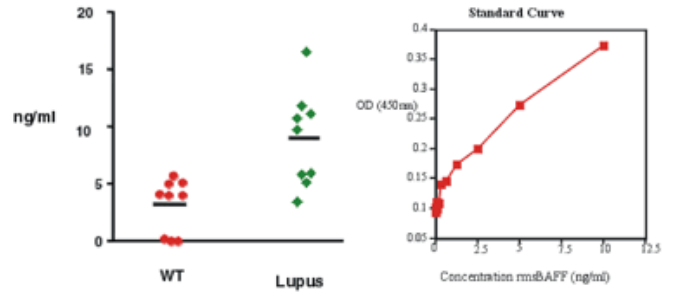


Figure: left) Serum levels of BAFF (mouse) in wild-type (red) and lupus-prone mice (green) measured with the matched pair MAbs 5A8 and 1C9. right) Standard curve: coated with 5A8 (5µg/ml) and detected with 1C9 (Biotin) followed by SAP-HRP.

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

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

Regulatory Status

RUO - Research Use Only

Product Details

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

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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