

BAFF (soluble) (mouse), (recombinant)

Citations: 3

[View Online »](#)

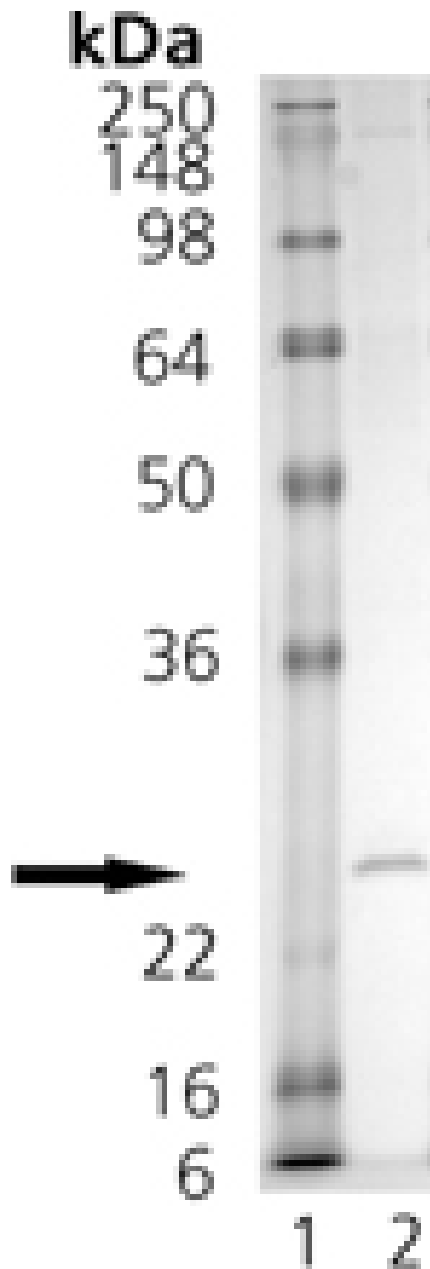
Ordering Information

[Order Online »](#)

| | |
|------------------|------|
| ALX-522-052-C010 | 10µg |
|------------------|------|

Manuals, SDS & CofA

[View Online »](#)



SDS-PAGE analysis: Lane 1: MW Marker, Lane 2: 1 μ g BAFF, stained with Imperial stain

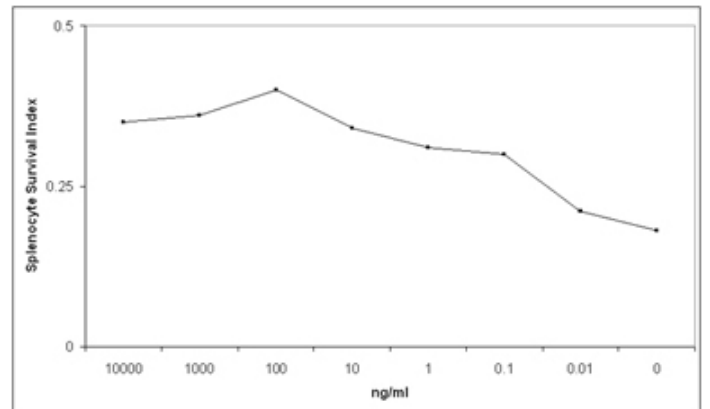


Figure: BAFF, Soluble (mouse) (rec.) mediates survival of freshly isolated splenocytes. Method: On day 0 splenocytes were isolated from a freshly collected C57Bl6 spleen. An aliquot of the splenocytes was analyzed on FACS and gated on the SSC-FSC panel. FACS settings were saved. The rest of the cells was put in culture with media alone or with increasing concentrations of BAFF (Prod. No. ALX-522-052) as indicated. After three days in culture, cells were harvested and analyzed on FACS with the saved setting. Splenocytes Survival Index (ratio % living/% dead cells) was calculated and plotted.

Handling & Storage

| | |
|-------------------|--|
| Use/Stability | Stable for at least 6 months after receipt when stored at -20°C. |
| Handling | Avoid freeze/thaw cycles. After reconstitution, prepare aliquots and store at -20°C. |
| 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 Notes | ELISA: binds to BAFF receptor at 1-100 ng/ml. |
| Concentration | 0.1mg/ml after reconstitution. |
| Endotoxin Content | <0.1EU/μg purified protein (LAL test; Associates of Cape Cod). |
| Formulation | Lyophilized. Contains PBS. |
| Gene/Protein Identifier | A000383 (UCSD Signaling Gateway ID) |
| MW | ~17kDa (SDS-PAGE). |
| Purity | ≥90% (SDS-PAGE) |
| Reconstitution | Reconstitute with 100μl sterile water. Further dilutions should be made with medium containing 5% fetal calf serum or other carrier protein. |
| Source | Produced in HEK 293 cells. The extracellular domain of mouse BAFF (aa 127-309) is fused at the N-terminus to a linker peptide (10 aa) and a FLAG [®] -tag. |
| Specificity | Binds to human (weak) and mouse BCMA, TACI and BAFF-R. |

Technical Info / Product Notes Historical data has shown that BAFF mediates splenocyte survival.

Notes



UniProt ID

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

FLAG is a registered trademark of Sigma-Aldrich Co.

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

ENZO LIFE SCIENCES & Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

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

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

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