

# BAFF (soluble) (mouse), (recombinant)

Citations: 3

[View Online »](#)

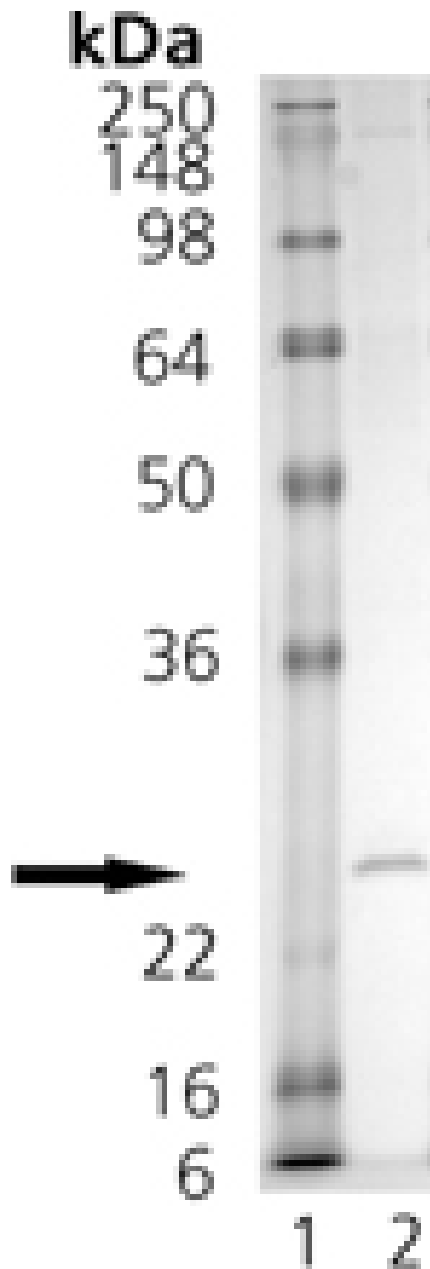
Ordering Information

[Order Online »](#)

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

Manuals, SDS & CofA

[View Online »](#)



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

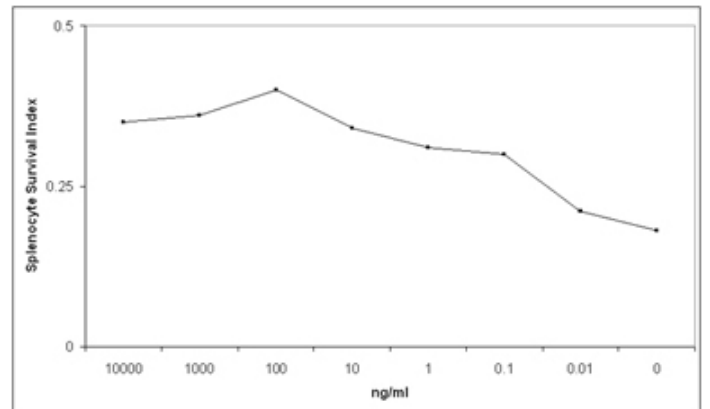


Figure: BAFF, Soluble (mouse) (rec.) mediates survival of freshly isolated splenocytes. Method: On day 0 splenocytes were isolated from a freshly collected C57Bl6 spleen. An aliquot of the splenocytes was analyzed on FACS and gated on the SSC-FSC panel. FACS settings were saved. The rest of the cells 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 <sup>®</sup> -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.

FLAG is a registered trademark of Sigma-Aldrich Co.

## UniProt ID

Q9WU72

Last modified: May 29, 2024



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-  
be@enzolifesciences.com](mailto:info-be@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)