

BAFF (soluble) (human), (recombinant)

Citations: 18

[View Online »](#)

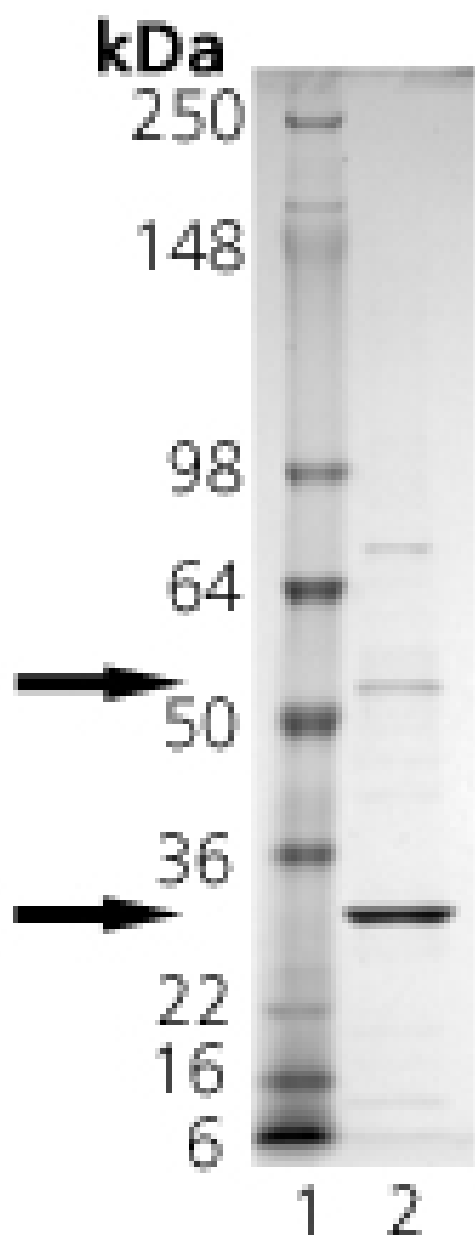
Ordering Information

[Order Online »](#)

ALX-522-025-C010	10µg
------------------	------

Manuals, SDS & CofA

[View Online »](#)



SDS-PAGE Analysis: Lane 1: MW Marker, Lane 2: 1 μ g BAFF. Bands indicated are immunoreactive with anti-FLAG antibody.

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	THANK, zTNF4, TALL-1, BlyS, TNFSF 13B/20, CD257
Application Notes	ELISA: binds to BAFF receptor at 2-100 ng/ml. Detection of receptors of human soluble BAFF on target cells.
Concentration	0.1mg/ml after reconstitution.
Endotoxin Content	<1EU/µg purified protein (LAL test; Associates of CAPE COD Inc.).
Formulation	Lyophilized. Contains PBS.
MW	~28kDa (SDS-PAGE).
Purity	≥95% (SDS-PAGE)
Reconstitution	Reconstitute with 100µl sterile water. Further dilutions should be made with medium containing 5% fetal calf serum.
Source	Produced in <i>E. coli</i> . The extracellular domain of human BAFF (aa 83-285) is fused at the N-terminus to a linker peptide (6 aa) and a FLAG [®] -tag.
Specificity	Binds to human and mouse (weak) BCMA, TACI and BAFF-R.

Technical Info / Product Notes

Historical data has shown that BAFF mediates splenocyte survival and induces monocyte survival.

FLAG is a registered trademark of Sigma-Aldrich Co.

UniProt ID

Q9Y275

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)