

BCMA (human):Fc (human), (recombinant)

Citations: 12

[View Online »](#)

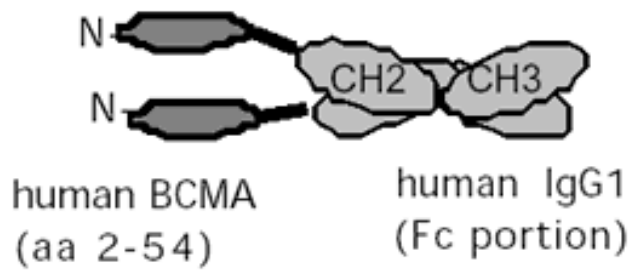
Ordering Information

[Order Online »](#)

ALX-522-026-C050	50µg
------------------	------

Manuals, SDS & CofA

[View Online »](#)



Schematic structure of human rhBCMA:Fc.

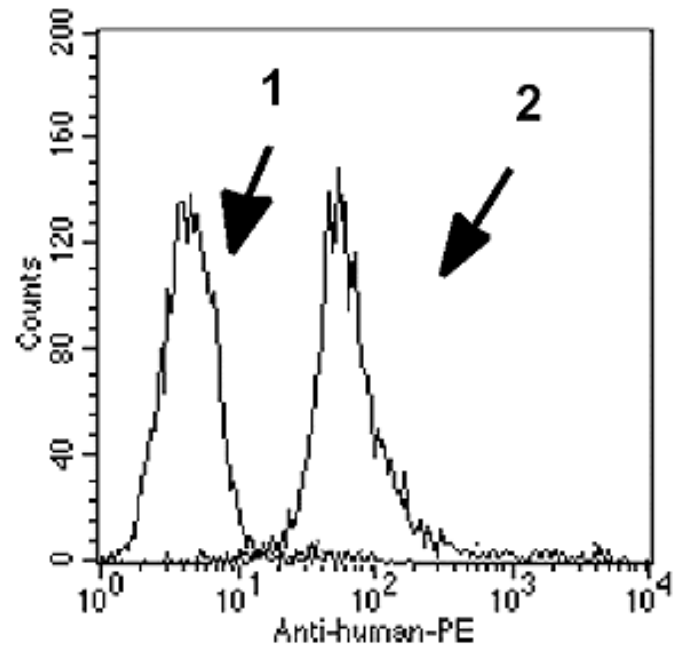


Figure: Flow cytometric profile of 293 cells stably transfected with a BAFF-expression plasmid (2). Control (1) are untransfected 293 cells. Optimal concentration range is 0.5-5 $\mu\text{g/ml}$. **Method:** 293 cells (5×10^5) expressing surface BAFF vector were incubated on ice for 30 min in 50 μl FACS buffer (PBS, 5% Fetal calf serum, 0.02% azide) containing 4 $\mu\text{g/ml}$ of BCMA:Fc. After washing in FACS buffer, PE-conjugated antibody to human IgG (SBA) was added and cells were analyzed by flow cytometry.

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	TNFRSF 17, B cell maturation protein, CD269
Application Notes	ELISA: binds to mouse BAFF). Flow Cytometry: Detection of membrane-bound mouse BAFF and human APRIL in combination with fluorescently labelled Ab to IgG1.
Concentration	1mg/ml after reconstitution.
Formulation	Lyophilized. Contains PBS.
MW	~40kDa (SDS-PAGE).
Purity	≥95% (SDS-PAGE)
Reconstitution	Reconstitute with 50µl of sterile water. Further dilutions should be made with medium containing 5% fetal calf serum or a carrier protein.
Source	Produced in HEK 293 cells. The extracellular domain of human BCMA (aa 2-54) is fused at the C-terminus to the Fc portion of human IgG1.
Specificity	Binds human and mouse BAFF and APRIL.
Technical Info / Product Notes	Historical data shows that it blocks the binding of BAFF and APRIL to their receptors BCMA and TACI, inhibiting BAFF- and APRIL-mediated B cell activation.

UniProt ID Q02223



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-be@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)