

BAFF-R (human):Fc (human), (recombinant)

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-522-060-C050	50µg
------------------	------

Manuals, SDS & CofA

[View Online »](#)

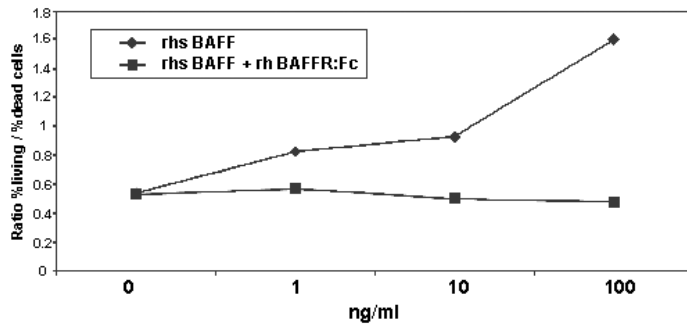


Figure: rhBAFF-R:Fc inhibits rhBAFF-mediated activation of freshly isolated mouse splenocytes. Concentration of rhBAFF-R:Fc required to achieve complete inhibition depends on the concentration of rhBAFF (Prod. No. [ALX-522-025](#)) used. A ratio (rhBAFF:rhBAFF-R:Fc) 1:50 and higher is recommended. **Method:** On day 0 splenocytes were isolated from a freshly collected C57B16 spleen. An aliquot of the splenocytes was analyzed on FACS, and the B-cells were gated on the SSC-FSC panel. FACS settings were saved. The rest of the cells was put in culture with media only, with the indicated concentration of rhBAFF (Prod. No. ALX-522-025) alone or in the presence of excess rhBAFF-R:Fc (+5µg/ml). After 3 days in culture

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 13C, BR3, BlySR3, CD268
Application Notes	ELISA: binds to BAFF. Detection of membrane-bound human and mouse BAFF in combination with PAb to human IgG1.
Concentration	1mg/ml (after reconstitution)
Endotoxin Content	<0.1EU/µg purified protein (LAL test; Bio Whittaker).
Formulation	Lyophilized. Contains PBS.
MW	~40kDa (SDS-PAGE)
Purity	≥95% (SDS-PAGE)
Reconstitution	Reconstitute with 50µl 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 BAFF-R (aa 2-71) is fused at the C-terminus to the Fc portion of human IgG1.
Specificity	Binds human and mouse BAFF.
Technical Info / Product Notes	Historical results have shown that BAFF-R inhibits rhsBAFF mediated splenocyte survival.
UniProt ID	Q96RJ3



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)