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

Citations: 1

View Online »

Ordering Information

Order Online »

ALX-522-060-C050

50µg

Manuals, SDS & CofA

View Online »

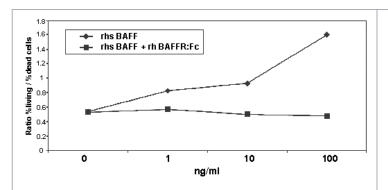


Figure: rhBAFF-R:Fc inhibits rhsBAFF-mediated activation of freshly isolated mouse splenocytes. Concentration of rhBAFF-R:Fc required to achieve complete inhibition depends on the concentration of rhsBAFF (Prod. No. ALX-522-025) used. A ratio (rhsBAFF:rhBAFF-R:Fc) 1:50 and higher is recommended.Method: 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 rhsBAFF (Prod. No. ALX-522-025) alone or in the presence of excess rhsBAFF-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

Last modified: May 29, 2024



info-

eu@enzolifesciences.com