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

Citations: 12

View Online »

Ordering Information

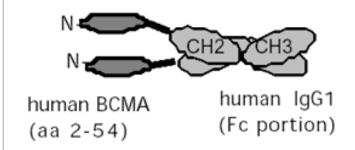
Order Online »

ALX-522-026-C050

50µg

Manuals, SDS & CofA

View Online »



Schematic structure of human rhBCMA:Fc.

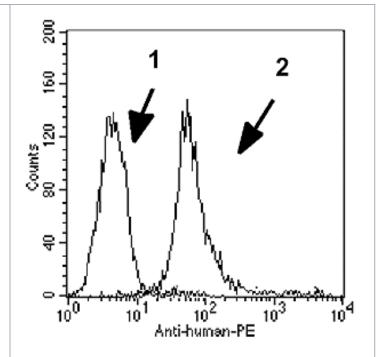


Figure: Flow cytometric profile of 293 cells stably transfected with a BAFF-expression plasmid (2). Control (1) are untransfected 293 cells. Optimal concentration range is 0.5-5 μ g/ml.**Method**: 293 cells (5 x 10⁵) expressing surface BAFF vector were incubated on ice for 30 min in 50 μ l FACS buffer (PBS, 5% Fetal calf serum, 0.02% azide) containing 4 μ 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.

SpecificityBinds human and mouse BAFF and APRIL.

Technical Info / Product NotesHistorical 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

Last modified: May 29, 2024



info-

eu@enzolifesciences.com