

XFD750 Anti-human CD23 Antibody *EBVCS-5*

Catalog Number: 102301A0, 102301A1

Unit Size: 100 tests, 500 tests

Product Details

Storage Conditions	2-8°C with minimized light exposure. Do not freeze.
Expiration Date	12 months upon receiving
Concentration	Lot specific (please consult certificate of analysis for given lot)
Formulation	Phosphate-buffered saline (PBS, pH 7.2), 0.09% sodium azide, 0.2% (w/v) BSA

Antibody Properties

Species Reactivity	Human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Isotype	Mouse IgG1, κ
Immunogen	CD23 (Fc ϵ RII, B6, BLAST-2, Leu-20, Low affinity IgE receptor)
Clone	EBVCS-5
Conjugate	AF750

Biological Properties

Appearance	Dark blue liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with AF750 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging
Recommended Dilutions	For flow cytometry applications, the suggested concentration is at 5 μ L/million cells in 100 μ L staining buffer. For the best performance of each application, the optimal concentration of this reagent needs to be carefully determined. <i>*The suggested working dilution is provided as a guide only. It is recommended that the users titrates the product for use in their tests using proper positive and negative controls.</i>

Spectral Properties

Conjugate	AF750
Excitation Wavelength	752 nm
Emission Wavelength	776 nm

Applications

EBVCS-5 is an anti-human antibody that forms an immune complex with the CD23 antigen. CD23 (sometimes referred to as FcεRII, BLAST-2 or B6) is a glycoprotein that is expressed on the surface of cells like platelets, granulocytes, T cells, dendritic cells and epithelial cells. CD23 is a component of important cellular pathways, for example, the cytokine-mediated signaling pathway and Notch signaling pathway. Furthermore, in certain organisms, it is a promoter of nitric-oxide synthase activity, acts to positively regulate killing of cells of other organism and positively regulates nitric-oxide synthase biosynthetic process. From a research standpoint, it is of biological interest due to its association with key macromolecules/ligands like CD11b, CD21, IgE and CD11c. CD23 is a fairly uncommon antibody target, with a little more than 5800 publications in the last decade. Even still, CD23 is typically used in flow cytometry applications as a phenotypic marker for differentiation of cell types, specifically in the study of . This antibody was purified through affinity chromatography and conjugated to XFD750 (ex/em = 752/776 nm). XFD750 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 750 (Alexa Fluor® is the trademark of Thermo Fisher).