

XFD750 Anti-human CD50 Antibody *MEM-04*

Catalog Number: 105001B0,

105001B1

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 IgG2b kappa
Immunogen	CD50 (ICAM-3)
Clone	MEM-04
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 uL/million cells in 100 uL 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

MEM-04 is an anti-human monoclonal antibody that targets the CD50 antigen. CD50 (alternatively called ICAM-3) is a 130 kD member of the Ig superfamily that is located on the surface of cells like macrophages. CD50 is associated with a variety of biologically interesting macromolecules/ligands, in particular, integrin α _d/β₂ and LFA-1. CD50 is a relatively rare antibody target, with fewer than 300 publications in the last decade. Even still, CD50 is typically used in flow cytometry applications as a phenotypic marker for differentiation of cell types, particularly in the study of costimulatory molecules and cell biology. 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).