

PE/XFD700 Anti-human CD50 Antibody *MEM-04*

Catalog Number: 105001Q0,
105001Q1, 105001Q2
Unit Size: 25 tests, 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	PE/AF700

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PE/AF700 under optimal conditions
Application	Flow Cytometry (FACS)
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	PE/AF700
-----------	----------

Excitation Wavelength 565 nm

Emission Wavelength 721 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 β 2 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 PE/XFD700 (ex/em = 566/721 nm). XFD700 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 700 (Alexa Fluor® is the trademark of Thermo Fisher). It is compatible with the 561 nm laser and 702/85 nm bandpass filter (for example, as in the Luminex Amnis FlowSight).