

PE/Cy7 Anti-human CD50 Antibody
MEM-04Catalog number: 10500100, 10500101, 10500102
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	0.1 mg/mL
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/Cy7

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PE/Cy7 under optimal conditions
Application	Flow Cytometry (FACS)

Spectral Properties

Conjugate	PE/Cy7
Excitation Wavelength	566 nm
Emission Wavelength	778 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/Cy7 (ex/em = 566/778 nm). It is compatible with the 561 nm laser and 780/60 nm bandpass filter (for example, as in the Agilent

Technologies NovoCyte Advanteon).