

PerCP Anti-human CD50 Antibody
MEM-04Catalog number: 105001V0, 105001V1, 105001V2
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	PerCP

Biological Properties

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

Spectral Properties

Conjugate	PerCP
Excitation Wavelength	477 nm
Emission Wavelength	678 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/ 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 PerCP (ex/em = 477/678 nm). It is compatible with the 488 nm laser and 693/37 nm bandpass filter (for example, as in the Miltenyi Biotec

MACSQuant Analyzer 10).