

**mFluor™ Green 620 Anti-human CD50
Antibody *MEM-04***Catalog number: 105000U0, 105000U1
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	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	mFluor™ Green 620

Biological Properties

Appearance	Purple liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with mFluor™ Green 620 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	mFluor™ Green 620
Excitation Wavelength	525 nm
Emission Wavelength	623 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 mFluor™ Green 620 (ex/em = 525/623 nm). It is compatible with the 532 nm laser and 611/31 nm bandpass filter (for example, as in the Luminex Amnis CellStream).