

**mFluor™ Violet 540 Anti-non-human
primates/ human CD48 Antibody
*MEM-102***Catalog number: 10480120, 10480121
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	Non-human primates, human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Isotype	Mouse IgG1
Immunogen	CD48 (Blast-1, Hulym3, BCM1)
Clone	MEM-102
Conjugate	mFluor™ Violet 540

Biological Properties

Appearance	Yellow liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with mFluor™ Violet 540 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	mFluor™ Violet 540
Excitation Wavelength	394 nm
Emission Wavelength	537 nm

Applications

MEM-102 is an anti-non-human primates/ human monoclonal antibody that targets the CD48 antigen. CD48 (sometimes referred to as Blast-1) is a 40 - 47 kD member of the CD2 family that is expressed on the surface of cells like T cells and dendritic cells. CD48 is associated with a variety

of biologically interesting macromolecules/ligands, for instance, CD229, fyn, CD244 and CD2. CD48 is a fairly uncommon antibody target, with a little more than 2800 publications in the last decade. Even still, CD48 is commonly used in flow cytometry applications as a phenotypic marker for differentiation of cell types, particularly in the study of immunology. This antibody was purified through affinity chromatography and conjugated to mFluor™ Violet 540 (ex/em = 394/537 nm). It is compatible with the 405 nm laser and 525/50 nm bandpass filter (for example, as in the BD FACSymphony™ A5).