

**iFluor™ 555 Anti-non-human primates/
human CD48 Antibody *MEM-102***Catalog number: 10480090, 10480091
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	iFluor™ 555

Biological Properties

Appearance	Red liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 555 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 555
Excitation Wavelength	557 nm
Emission Wavelength	570 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 iFluor™ 555 (ex/em = 557/570 nm). It is compatible with the 561 nm laser and 586/14 nm bandpass filter (for example, as in the BD LSRFortessa™ X-20).