

**mFluor™ Blue 570 Anti-human CD68  
Antibody \*Y1/82A\***Catalog number: 106800T0, 106800T1  
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 IgG2a
Immunogen	CD68 (Macrosialin)
Clone	Y1/82A
Conjugate	mFluor™ Blue 570

**Biological Properties**

Appearance	Red liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with mFluor™ Blue 570 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

**Spectral Properties**

Conjugate	mFluor™ Blue 570
Excitation Wavelength	553 nm
Emission Wavelength	565 nm

**Applications**

The Y1/82A monoclonal antibody recognizes human CD68, a 110 kD glycoprotein typically located on the surface of T cells, Iak cells, dendritic cells, B cells and neutrophils. CD68 is associated with a variety of biologically interesting macromolecules/ligands, namely, LDL. CD68 is a very popular antibody target, with over 21000 publications in the last decade. CD68 is frequently used in flow cytometry applications as a phenotypic

marker for differentiation of cell types, particularly in the study of neuroscience cell markers, immunology and neuroscience. This antibody was purified through affinity chromatography and conjugated to mFluor™ Blue 570 (ex/em = 553/565 nm). It is compatible with the 561 nm laser and 582/15 nm bandpass filter (for example, as in the BD FACSAria™ III).