

TRITC Anti-human CD68 Antibody *Y1/82A*Catalog number: 106801J0, 106801J1
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	TRITC

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with TRITC under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	TRITC
Excitation Wavelength	544 nm
Emission Wavelength	570 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 TRITC (ex/em = 544/570 nm). It is compatible with the 561 nm laser and 586/20 nm bandpass filter (for example, as in the Agilent Technologies NovoCyte).