

iFluor™ 790 Anti-human CD68 Antibody
Y1/82ACatalog number: 106800M0, 106800M1
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	iFluor™ 790

Biological Properties

Appearance	Green liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 790 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 790
Excitation Wavelength	787 nm
Emission Wavelength	812 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 iFluor™ 790 (ex/em = 787/812 nm).