

iFluor™ 350 Anti-human CD64 Antibody
10.1Catalog number: 10640010, 10640011
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 IgG1
Immunogen	CD64 (FcR I)
Clone	10.1
Conjugate	iFluor™ 350

Biological Properties

Appearance	Off-white liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 350 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 350
Excitation Wavelength	345 nm
Emission Wavelength	450 nm

Applications

The 10.1 monoclonal antibody binds with human CD64, a 72 kD single-pass type I membrane protein typically expressed on the surface of granulocytes, monocytes and dendritic cells. In some organisms, CD64 enhances protein tyrosine kinase activity. Moreover, it is a member of vital cellular pathways, for example, the interferon-gamma-mediated signaling pathway and Fc-gamma receptor signaling pathway involved in

phagocytosis. From a research standpoint, it is of biological interest due to its association with important macromolecules/ligands such as . CD64 is a fairly uncommon antibody target, with a little more than 4000 publications in the last decade. Even still, CD64 is often used in flow cytometry applications as a phenotypic marker for differentiation of cell types, particularly in the study of immunology and innate immunity. This antibody was purified through affinity chromatography and conjugated to iFluor™ 350 (ex/em = 345/450 nm). It is compatible with the 355 nm laser and 447/60 nm bandpass filter (for example, as in the Bio-Rad ZE5 Cell Analyzer).