

iFluor™ 546 Anti-human CD11b Antibody
ICRF44Catalog number: 10112080, 10112081
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, κ
Immunogen	CD11b (CR3, Mac-1, Mo1, ITGAM, Integrin alpha-M)
Clone	ICRF44
Conjugate	iFluor™ 546

Biological Properties

Appearance	Red liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 546 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 546
Excitation Wavelength	541 nm
Emission Wavelength	557 nm

Applications

ICRF44 is an anti-human monoclonal antibody that forms an immune complex with the CD11b antigen. CD11b (also known as Mo1 or Mac-1) is a 165 - 170 kD transmembrane glycoprotein that is expressed on the surface of cells such as dendritic cells, macrophages, NK cells and T cells. In some organisms, CD11b is involved in the positive regulation of neuron death, is a promoter of hippocampal neuron apoptotic process and is a

positive regulator of protein targeting to membrane. Furthermore, it has been thought to be involved with vital biological processes such as cell adhesion, specifically cell-cell adhesion via plasma-membrane adhesion molecules. CD11b is a member of vital cellular pathways, for example, the integrin-mediated signaling pathway, apoptotic signaling pathway and toll-like receptor 4 signaling pathway. From a research standpoint, it is of biological interest due to its association with essential macromolecules/ligands like iC3b, ICAM-1, Factor X and 2. CD11b is a very popular antibody target, with over 45000 publications in the last decade. CD11b is frequently used in flow cytometry applications as a phenotypic marker for differentiation of cell types, especially in the study of innate immunity and immunology. This antibody was purified through affinity chromatography and conjugated to iFluor™ 546 (ex/em = 541/557 nm). It is compatible with the 532 nm laser and 575/36 nm bandpass filter (for example, as in the Thermo Fisher Attune NxT).