

**mFluor™ Red 780 Anti-human CD11b  
Antibody \*ICRF44\***Catalog number: 101120W0, 101120W1  
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          | mFluor™ Red 780                                  |

**Biological Properties**

|             |                                                                                                                |
|-------------|----------------------------------------------------------------------------------------------------------------|
| Appearance  | Dark blue liquid                                                                                               |
| Preparation | Antibody purified by affinity chromatography and then conjugated with mFluor™ Red 780 under optimal conditions |
| Application | Flow Cytometry (FACS), Fluorescence Imaging                                                                    |

**Spectral Properties**

|                       |                 |
|-----------------------|-----------------|
| Conjugate             | mFluor™ Red 780 |
| Excitation Wavelength | 629 nm          |
| Emission Wavelength   | 767 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 mFluor™ Red 780 (ex/em = 629/767 nm). It is compatible with the 633 nm laser and 783/56 nm bandpass filter (for example, as in the BD FACSVerser™).