

**mFluor™ Red 700 Anti-human CD11c  
Antibody \*3.9\***Catalog number: 101130V0, 101130V1  
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          | CD11c (Integrin alpha-X, CR4, p150, ITGAX) |
| Clone              | 3.9                                        |
| Conjugate          | mFluor™ Red 700                            |

**Biological Properties**

|             |                                                                                                                |
|-------------|----------------------------------------------------------------------------------------------------------------|
| Appearance  | Dark blue liquid                                                                                               |
| Preparation | Antibody purified by affinity chromatography and then conjugated with mFluor™ Red 700 under optimal conditions |
| Application | Flow Cytometry (FACS), Fluorescence Imaging                                                                    |

**Spectral Properties**

|                       |                 |
|-----------------------|-----------------|
| Conjugate             | mFluor™ Red 700 |
| Excitation Wavelength | 680 nm          |
| Emission Wavelength   | 695 nm          |

**Applications**

3.9 is an anti-human monoclonal antibody that forms an immune complex with the CD11c antigen. CD11c (sometimes called CR4 or ITGAX) is a 145 - 150 kD single-pass type I membrane protein that is expressed on the surface of cells such as macrophages, B cells and dendritic cells. In some organisms, CD11c promotes angiogenesis, positively regulates myelination and is a promoter of endothelial tube morphogenesis.

Moreover, it is a component of essential cellular pathways, in particular, the cytokine-mediated signaling pathway and integrin-mediated signaling pathway. From a research standpoint, it is of biological interest due to its association with vital macromolecules/ligands like ICAM-1 and 4 and fibrinogen. CD11c is a very popular antibody target, with over 26000 publications in the last decade. CD11c has been widely used in costimulatory molecules research, commonly serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to mFluor™ Red 700 (ex/em = 680/695 nm). It is compatible with the 642 nm laser and 664/20 nm bandpass filter (for example, as in the Luminex Guava easyCyte).