

**PacBlue Anti-human CD11c Antibody \*3.9\***Catalog number: 101131J0, 101131J1  
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, $\kappa$
Immunogen	CD11c (Integrin alpha-X, CR4, p150, ITGAX)
Clone	3.9
Conjugate	PacBlue

**Biological Properties**

Appearance	Light yellow liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with PacBlue under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

**Spectral Properties**

Conjugate	PacBlue
Excitation Wavelength	404 nm
Emission Wavelength	455 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 PacBlue (ex/em = 404/455 nm). It is compatible with the 405 nm laser and 450/45 nm bandpass filter (for example, as in the Luminex Guava easyCyte).