

PE/iFluor™ 700 Anti-human CD11c Antibody
3.9Catalog number: 101131X0, 101131X1, 101131X2
Unit size: 25 tests, 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	PE/iFluor™ 700

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PE/iFluor™ 700 under optimal conditions
Application	Flow Cytometry (FACS)

Spectral Properties

Conjugate	PE/iFluor™ 700
Excitation Wavelength	566 nm
Emission Wavelength	708 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 PE/iFluor™ 700 (ex/em = 566/708 nm). It is compatible with the 561 nm laser and 695/40 nm bandpass filter (for example, as in the Agilent Technologies NovoCyte).