

PE/Cy7 Anti-human CD11b Antibody
HI11bCatalog number: 101111N0, 101111N1, 101111N2
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 IgG2b
Immunogen	CD11b (CR3, Mac-1, Mo1, ITGAM, Integrin alpha-M)
Clone	HI11b
Conjugate	PE/Cy7

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PE/Cy7 under optimal conditions
Application	Flow Cytometry (FACS)

Spectral Properties

Conjugate	PE/Cy7
Excitation Wavelength	566 nm
Emission Wavelength	778 nm

Applications

HI11b is an anti-human monoclonal antibody that recognizes the CD11b antigen. CD11b (sometimes referred to as Integrin α M chain, C3bIR, CR3A or CR3) is a 165 - 170 kD glycoprotein that is expressed on the surface of cells such as B cells, granulocytes, NK cells, macrophages and dendritic cells. CD11b is involved with important cellular pathways, in particular, the cytokine-mediated signaling pathway, toll-like receptor 4 signaling pathway and apoptotic signaling pathway. Additionally, in many organisms, it is a promoter of superoxide anion generation, positively regulates microglial cell mediated cytotoxicity and is a positive regulator of prostaglandin-E synthase activity. CD11b has been associated with critical biological processes like cell adhesion, specifically cell-cell adhesion via plasma-membrane adhesion molecules, and is associated with a

variety of biologically interesting macromolecules/ligands, for example, Factor X, iC3b and ICAM-1. CD11b is a very popular antibody target, with over 45000 publications in the last decade. CD11b has been widely used in innate immunity, neuroscience cell markers and cell biology research, often serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to PE/Cy7 (ex/em = 566/778 nm). It is compatible with the 561 nm laser and 780/60 nm bandpass filter (for example, as in the BD FACSCelesta™).