

**iFluor™ 405 Anti-human CD11b Antibody**  
**\*HI11b\***Catalog number: 10111020, 10111021  
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 IgG2b
Immunogen	CD11b (CR3, Mac-1, Mo1, ITGAM, Integrin alpha-M)
Clone	HI11b
Conjugate	iFluor™ 405

**Biological Properties**

---

Appearance	Light yellow liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 405 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

**Spectral Properties**

---

Conjugate	iFluor™ 405
Excitation Wavelength	403 nm
Emission Wavelength	427 nm

**Applications**

---

HI11b is an anti-human monoclonal antibody that recognizes the CD11b antigen. CD11b (sometimes referred to as Integrin  $\alpha$ 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 iFluor™ 405 (ex/em = 403/427 nm). It is compatible with the 405 nm laser and 445/45 nm bandpass filter (for example, as in the Agilent Technologies NovoCyte Quanteon).