

**PerCP/Cy5.5 Anti-human CD11a Antibody**  
**\*HI111\***Catalog number: 101101U0, 101101U1, 101101U2  
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
Immunogen	CD11a (LFA-1A, Integrin $\alpha$ L, ITGAL)
Clone	HI111
Conjugate	PerCP/Cy5.5

**Biological Properties**

Preparation	Antibody purified by affinity chromatography and then conjugated with PerCP/Cy5.5 under optimal conditions
Application	Flow Cytometry (FACS)

**Spectral Properties**

Conjugate	PerCP/Cy5.5
Excitation Wavelength	489 nm
Emission Wavelength	679 nm

**Applications**

HI111 is an anti-human monoclonal antibody that forms an immune complex with the CD11a antigen. CD11a (also known as Integrin  $\alpha$ -L or LFA-1 $\alpha$  chain) is a 170 - 180 kD transmembrane protein that is expressed on the surface of cells such as granulocytes and macrophages. CD11a has been associated with critical biological processes like cell-cell adhesion, especially leukocyte cell-cell adhesion. Moreover, it is a member of essential cellular pathways, for instance, the integrin-mediated signaling pathway. From a research standpoint, it is of biological interest due to its association with vital macromolecules/ligands such as CD18 and ICAM-1, 2, 3 and 4. CD11a is a fairly uncommon antibody target, with a little

more than 3700 publications in the last decade. Even still, CD11a has a variety of applications in neuroscience and innate immunity research, typically serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to PerCP/Cy5.5 (ex/em = 489/679 nm). It is compatible with the 488 nm laser and 700/54 nm bandpass filter (for example, as in the BD FACSMelody™).