

**PerCP/Cy5.5 Anti-human CD82 Antibody**  
**\*C33\***Catalog number: 108201W0, 108201W1, 108201W2  
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 IgG2a
Immunogen	CD82 (KAI1, C33, R2, 4F9, IA4)
Clone	C33
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**

---

The C33 monoclonal antibody binds with human CD82, a 45 - 90 kD member of the 4-span transmembrane protein superfamily often located on the surface of epithelial cells, fibroblasts, endothelial cells and lymphocytes. CD82 is associated with a variety of biologically interesting macromolecules/ligands, in particular, MHC-II. CD82 is a fairly uncommon antibody target, with a little more than 1600 publications in the last decade. Even still, CD82 is essential for immunology 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 PerCP/Cy5.5 (ex/em = 489/679 nm). It is

compatible with the 488 nm laser and 690/50 nm bandpass filter (for example, as in the Beckman Coulter DxFLEx).