

PerCP Anti-human CD82 Antibody *C33*

Catalog number: 108201V0, 108201V1, 108201V2

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

Biological Properties

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

Spectral Properties

Conjugate	PerCP
Excitation Wavelength	477 nm
Emission Wavelength	678 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 (ex/em = 477/678 nm). It is compatible with the 488 nm laser and 695/40 nm bandpass filter (for example, as in the BD FACSAria™ Fusion).