

iFluor™ 568 Anti-human CD82 Antibody
C33Catalog number: 108200B0, 108200B1
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 IgG2a
Immunogen	CD82 (KAI1, C33, R2, 4F9, IA4)
Clone	C33
Conjugate	iFluor™ 568

Biological Properties

Appearance	Purple liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 568 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 568
Excitation Wavelength	568 nm
Emission Wavelength	587 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 iFluor™ 568 (ex/em = 568/587 nm). It is compatible with the 561 nm laser and 585/29 nm bandpass filter (for example, as in the BD FACSJazz™).