

iFluor™ 670 Anti-human CD100 Antibody
133-1C6Catalog number: 110000H0, 110000H1
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 IgM
Immunogen	CD100 (SEMA4D)
Clone	133-1C6
Conjugate	iFluor™ 670

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 670 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 670
Excitation Wavelength	671 nm
Emission Wavelength	682 nm

Applications

The 133-1C6 monoclonal antibody recognizes human CD100, a 150 kD member of the Class IV semaphorin family frequently expressed on the surface of B cells, T cells, oligodendrocytes and leukocytes. CD100 is associated with a variety of biologically interesting macromolecules/ligands, for instance, CD45. CD100 is a relatively rare antibody target, with fewer than 400 publications in the last decade. Even still, CD100 is frequently used in flow cytometry applications as a phenotypic marker for differentiation of cell types, particularly in the study of immunology. This antibody was purified through affinity chromatography and conjugated to iFluor™ 670 (ex/em = 671/682 nm). It is compatible

with the 642 nm laser and 664/20 nm bandpass filter (for example, as in the Luminex Guava easyCyte).