

**iFluor™ 670 Anti-non-human primates/
human CD170 Antibody *1A5***Catalog number: 117000H0, 117000H1
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	Non-human primates, human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Isotype	Mouse IgG1
Immunogen	CD170 (SIGLEC5, OB-BP2)
Clone	1A5
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

1A5 is an anti-non-human primates/ human monoclonal antibody that recognizes the CD170 antigen. CD170 (alternatively called OB-BP2 or SIGLEC5) is a 140 kD transmembrane protein that is expressed on the surface of cells such as dendritic cells, granulocytes, B cells and macrophages. CD170 is associated with a variety of biologically interesting macromolecules/ligands, namely, sialylated glycans. CD170 is a relatively rare antibody target, with fewer than 70 publications in the last decade. Even still, CD170 is commonly 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 702/87 nm bandpass filter (for example, as in the Luminex Amnis CellStream).