

**iFluor™ 555 Anti-non-human primates/  
human CD170 Antibody \*1A5\***Catalog number: 11700090, 11700091  
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™ 555               |

**Biological Properties**

|             |                                                                                                            |
|-------------|------------------------------------------------------------------------------------------------------------|
| Appearance  | Red liquid                                                                                                 |
| Preparation | Antibody purified by affinity chromatography and then conjugated with iFluor™ 555 under optimal conditions |
| Application | Flow Cytometry (FACS), Fluorescence Imaging                                                                |

**Spectral Properties**

|                       |             |
|-----------------------|-------------|
| Conjugate             | iFluor™ 555 |
| Excitation Wavelength | 557 nm      |
| Emission Wavelength   | 570 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™ 555 (ex/em = 557/570 nm). It is compatible with the 561 nm laser and 583/24 nm bandpass filter (for example, as in the Luminex Amnis CellStream).