

**iFluor™ 594 Anti-human CD99 Antibody**  
**\*HI156\***Catalog number: 109900C0, 109900C1  
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          | CD99 (MIC2, E2 antigen, HBA71, MSK5X) |
| Clone              | HI156                                 |
| Conjugate          | iFluor™ 594                           |

**Biological Properties**

|             |                                                                                                            |
|-------------|------------------------------------------------------------------------------------------------------------|
| Appearance  | Purple liquid                                                                                              |
| Preparation | Antibody purified by affinity chromatography and then conjugated with iFluor™ 594 under optimal conditions |
| Application | Flow Cytometry (FACS), Fluorescence Imaging                                                                |

**Spectral Properties**

|                       |             |
|-----------------------|-------------|
| Conjugate             | iFluor™ 594 |
| Excitation Wavelength | 588 nm      |
| Emission Wavelength   | 604 nm      |

**Applications**

The HI156 monoclonal antibody reacts with human CD99, a glycoprotein frequently found on the surface of endothelial cells, epithelial cells, thymocytes, natural killer cells and eosinophils. In certain organisms, CD99 positively regulates neutrophil extravasation, and is associated with a variety of biologically interesting macromolecules/ligands, for instance, . CD99 is a fairly uncommon antibody target, with a little more than

3000 publications in the last decade. Even still, CD99 has been widely used in immunology research, typically serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to iFluor™ 594 (ex/em = 588/604 nm). It is compatible with the 592 nm laser and 610/30 nm bandpass filter (for example, as in the Luminex Amnis ImageStream).