

**APC/iFluor™ 700 Anti-human CD99  
Antibody \*HI156\***Catalog number: 109901F0, 109901F1, 109901F2  
Unit size: 25 tests, 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	APC/iFluor™ 700

**Biological Properties**

Preparation	Antibody purified by affinity chromatography and then conjugated with APC/iFluor™ 700 under optimal conditions
Application	Flow Cytometry (FACS)

**Spectral Properties**

Conjugate	APC/iFluor™ 700
Excitation Wavelength	685 nm
Emission Wavelength	710 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

APC/iFluor™ 700 (ex/em = 685/710 nm).