

**APC Anti-human CD99 Antibody \*HIT4\***

Catalog number: 109911C0, 109911C1, 109911C2

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 IgM
Immunogen	CD99 (MIC2, E2 antigen, HBA71, MSK5X)
Clone	HIT4
Conjugate	APC

**Biological Properties**

---

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

**Spectral Properties**

---

Conjugate	APC
Excitation Wavelength	651 nm
Emission Wavelength	660 nm

**Applications**

---

The HIT4 monoclonal antibody reacts with human CD99, a transmembrane glycoprotein commonly located on the surface of epithelial cells, t cells, lymphocytes, natural killer cells and monocytes. In some 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 is often used in flow cytometry applications as a phenotypic marker for differentiation of cell types, especially in the study of immunology. This antibody was purified through affinity chromatography and conjugated to APC (ex/em = 651/660 nm). It is compatible with the 642 nm laser and 702/85 nm bandpass filter (for example, as in the Luminex Amnis FlowSight).