

**APC/iFluor™ 750 Anti-human CD99
Antibody *HIT4***Catalog number: 109911G0, 109911G1, 109911G2
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/iFluor™ 750

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with APC/iFluor™ 750 under optimal conditions
Application	Flow Cytometry (FACS)

Spectral Properties

Conjugate	APC/iFluor™ 750
Excitation Wavelength	754 nm
Emission Wavelength	793 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/iFluor™

750 (ex/em = 754/793 nm).