

iFluor™ 405 Anti-human CD99 Antibody
HIT4Catalog number: 10991020, 10991021
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 IgM
Immunogen	CD99 (MIC2, E2 antigen, HBA71, MSK5X)
Clone	HIT4
Conjugate	iFluor™ 405

Biological Properties

Appearance	Light yellow liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 405 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 405
Excitation Wavelength	403 nm
Emission Wavelength	427 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 iFluor™ 405 (ex/em = 403/427 nm). It is compatible with the 405 nm laser and 457/45 nm bandpass filter (for example, as in the Luminex Amnis FlowSight).