

PE/XFD610 Anti-human CD99 Antibody *HI156*

Catalog Number: 109901O0,
109901O1, 109901O2
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	Lot specific (please consult certificate of analysis for given lot)
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	PE/AF610

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PE/AF610 under optimal conditions
Application	Flow Cytometry (FACS)
Recommended Dilutions	For flow cytometry applications, the suggested concentration is at 5 uL/million cells in 100 uL staining buffer. For the best performance of each application, the optimal concentration of this reagent needs to be carefully determined.
	<i>*The suggested working dilution is provided as a guide only. It is recommended that the users titrates the product for use in their tests using proper positive and negative controls.</i>

Spectral Properties

Conjugate	PE/AF610
-----------	----------

Excitation Wavelength 565 nm

Emission Wavelength 627 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 PE/XFD610 (ex/em = 567/627 nm). XFD610 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 610 (Alexa Fluor® is the trademark of Thermo Fisher). It is compatible with the 561 nm laser and 610/20 nm bandpass filter (for example, as in the BD FACSAria™ Fusion).