

PE/XFD700 Anti-human CD43 Antibody *HI161*

Catalog Number: 104311P0,
104311P1, 104311P2
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 IgG1
Immunogen	CD43 (Leukocyte Sialoglycoprotein, Leukosialin, Galactoglycoprotein, SPN)
Clone	HI161
Conjugate	PE/AF700

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PE/AF700 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/AF700
-----------	----------

Excitation Wavelength 565 nm

Emission Wavelength 721 nm

Applications

The HI161 monoclonal antibody reacts with human CD43, a 95 - 135 kD transmembrane protein commonly located on the surface of plasma cells, thymocytes, neutrophils, myelomas and T cells. In some organisms, CD43 plays a role in the downregulation of cell adhesion, promotes tumor necrosis factor biosynthetic process and acts to negatively regulate T cell proliferation. From a research standpoint, it is of biological interest due to its association with vital macromolecules/ligands such as EZR. CD43 is a fairly uncommon antibody target, with a little more than 5000 publications in the last decade. Even still, CD43 has been widely used in immunology research, frequently serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to PE/XFD700 (ex/em = 566/721 nm). XFD700 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 700 (Alexa Fluor® is the trademark of Thermo Fisher). It is compatible with the 561 nm laser and 702/85 nm bandpass filter (for example, as in the Luminex Amnis FlowSight).