

XFD700 Anti-human CD71 Antibody *HI160*

Catalog Number: 107101A0, 107101A1

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	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 IgG2b
Immunogen	CD71 (T9, Transferrin receptor, TFRC)
Clone	HI160
Conjugate	AF700

Biological Properties

Appearance	Dark blue liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with AF700 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging
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	AF700
Excitation Wavelength	696 nm
Emission Wavelength	719 nm

Applications

HI160 is an anti-human monoclonal antibody that is specific for the CD71 antigen. CD71 (also known as Transferrin receptor protein 1) is a 95 kD member of the Transferrin receptor family that is located on the surface of cells such as endothelial cells and stem cells. In certain organisms, CD71 is a positive regulator of isotype switching, is involved in the positive regulation of T cell proliferation and acts to positively regulate B cell proliferation. From a research standpoint, it is of biological interest due to its association with important macromolecules/ligands such as Transferrin. CD71 is a fairly uncommon antibody target, with a little more than 4000 publications in the last decade. Even still, CD71 is commonly 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 XFD700 (ex/em = 696/719 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).