

PE/iFluor™ 700 Anti-human CD71 Antibody
HI160Catalog number: 107101X0, 107101X1, 107101X2
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 IgG2b
Immunogen	CD71 (T9, Transferrin receptor, TFRC)
Clone	HI160
Conjugate	PE/iFluor™ 700

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PE/iFluor™ 700 under optimal conditions
Application	Flow Cytometry (FACS)

Spectral Properties

Conjugate	PE/iFluor™ 700
Excitation Wavelength	566 nm
Emission Wavelength	708 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 PE/iFluor™ 700 (ex/em = 566/708 nm). It is compatible with the 561 nm laser and 695/40 nm bandpass filter (for example, as in the Agilent Technologies NovoCyte Advanteon).