

**iFluor™ 450 Anti-human CD71 Antibody**  
**\*HI160\***Catalog number: 10710040, 10710041  
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 IgG2b
Immunogen	CD71 (T9, Transferrin receptor, TFRC)
Clone	HI160
Conjugate	iFluor™ 450

**Biological Properties**

Appearance	Brown liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 450 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

**Spectral Properties**

Conjugate	iFluor™ 450
Excitation Wavelength	451 nm
Emission Wavelength	502 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 iFluor™ 450 (ex/em = 451/502 nm). It is compatible with the 445 nm laser and 510/80 nm bandpass filter (for example, as in the BD FACSAria™ III).