

iFluor™ 750 Anti-human CD71 Antibody
HI166Catalog number: 107110L0, 107110L1
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 IgG1
Immunogen	CD71 (T9, Transferrin receptor, TFRC)
Clone	HI166
Conjugate	iFluor™ 750

Biological Properties

Appearance	Dark blue liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 750 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 750
Excitation Wavelength	757 nm
Emission Wavelength	779 nm

Applications

HI166 is an anti-human monoclonal antibody that is specific for the CD71 antigen. CD71 (sometimes called Transferrin receptor or TFRC) 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 some organisms, CD71 is an enhancer of bone resorption, plays a role in the upregulation of B cell proliferation and is an enhancer of T cell proliferation. From a

research standpoint, it is of biological interest due to its association with essential 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 essential for immunology research, commonly serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to iFluor™ 750 (ex/em = 757/779 nm).