

**APC/iFluor™ 750 Anti-human CD71  
Antibody \*HI166\***Catalog number: 107111G0, 107111G1, 107111G2  
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 IgG1                            |
| Immunogen          | CD71 (T9, Transferrin receptor, TFRC) |
| Clone              | HI166                                 |
| Conjugate          | APC/iFluor™ 750                       |

**Biological Properties**

|             |                                                                                                                |
|-------------|----------------------------------------------------------------------------------------------------------------|
| Preparation | Antibody purified by affinity chromatography and then conjugated with APC/iFluor™ 750 under optimal conditions |
| Application | Flow Cytometry (FACS)                                                                                          |

**Spectral Properties**

|                       |                 |
|-----------------------|-----------------|
| Conjugate             | APC/iFluor™ 750 |
| Excitation Wavelength | 754 nm          |
| Emission Wavelength   | 793 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 APC/iFluor™ 750 (ex/em = 754/793 nm).