

**iFluor™ 568 Anti-human CD71 Antibody**  
**\*OKT-9\***Catalog number: 107130B0, 107130B1  
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, κ                         |
| Immunogen          | CD71 (T9, Transferrin receptor, TFRC) |
| Clone              | OKT-9                                 |
| Conjugate          | iFluor™ 568                           |

**Biological Properties**

|             |                                                                                                            |
|-------------|------------------------------------------------------------------------------------------------------------|
| Appearance  | Purple liquid                                                                                              |
| Preparation | Antibody purified by affinity chromatography and then conjugated with iFluor™ 568 under optimal conditions |
| Application | Flow Cytometry (FACS), Fluorescence Imaging                                                                |

**Spectral Properties**

|                       |             |
|-----------------------|-------------|
| Conjugate             | iFluor™ 568 |
| Excitation Wavelength | 568 nm      |
| Emission Wavelength   | 587 nm      |

**Applications**

The OKT-9 monoclonal antibody binds with human CD71, a 95 kD glycoprotein commonly expressed on the surface of reticulocytes, proliferating cells, reticulocytes and erythroid precursors. In many organisms, CD71 acts to positively regulate B cell proliferation, is a positive regulator of bone resorption and is an enhancer of T cell proliferation. From a research standpoint, it is of biological interest due to its association with vital

macromolecules/ligands like Transferrin. CD71 is a fairly uncommon antibody target, with a little more than 4300 publications in the last decade. Even still, CD71 is frequently 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™ 568 (ex/em = 568/587 nm). It is compatible with the 561 nm laser and 586/20 nm bandpass filter (for example, as in the Agilent Technologies NovoCyte).