

**iFluor™ 555 Anti-pig/ horse/ human/ non-human primates CD90 Antibody \*5E10\***Catalog number: 10900090, 10900091  
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 | Pig, horse, human, non-human primates |
| Class              | Primary                               |
| Clonality          | Monoclonal                            |
| Host               | Mouse                                 |
| Isotype            | Mouse IgG1 kappa                      |
| Immunogen          | CD90 (Thy-1)                          |
| Clone              | 5E10                                  |
| Conjugate          | iFluor™ 555                           |

**Biological Properties**

---

|             |  |
|-------------|--|
| Appearance  | Red liquid   |
| Preparation | Antibody purified by affinity chromatography and then conjugated with iFluor™ 555 under optimal conditions |
| Application | Flow Cytometry (FACS), Fluorescence Imaging  |

**Spectral Properties**

---

|                       |             |
|-----------------------|-------------|
| Conjugate             | iFluor™ 555 |
| Excitation Wavelength | 557 nm      |
| Emission Wavelength   | 570 nm      |

**Applications**

---

5E10 is an anti-pig/ horse/ human/ non-human primates monoclonal antibody that forms an immune complex with the CD90 antigen. CD90 (alternatively called Thy-1) is a 25 - 35 kD member of the Ig superfamily that is found on the surface of cells such as endothelial cells and stem cells. CD90 is associated with a variety of biologically interesting macromolecules/ligands, namely, fyn and p100. CD90 is a moderately popular

antibody target, with over 15000 publications in the last decade. CD90 is vital to immunology research, often serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to iFluor™ 555 (ex/em = 557/570 nm). It is compatible with the 561 nm laser and 582/15 nm bandpass filter (for example, as in the BD FACSMelody™).