

iFluor™ 350 Anti-pig/ horse/ human/ non-human primates CD90 Antibody *5E10*Catalog number: 10900010, 10900011
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™ 350 |

Biological Properties

| | |
|-------------|--|
| Appearance | Off-white liquid |
| Preparation | Antibody purified by affinity chromatography and then conjugated with iFluor™ 350 under optimal conditions |
| Application | Flow Cytometry (FACS), Fluorescence Imaging |

Spectral Properties

| | |
|-----------------------|-------------|
| Conjugate | iFluor™ 350 |
| Excitation Wavelength | 345 nm |
| Emission Wavelength | 450 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™ 350 (ex/em = 345/450 nm). It is compatible with the 355 nm laser and 447/60 nm bandpass filter (for example, as in the Bio-Rad ZE5 Cell Analyzer).