

# XFD700 Anti-pig/ horse/ human/ non-human primates CD90 Antibody \*5E10\*

Catalog Number: 109001A0, 109001A1

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	Lot specific (please consult certificate of analysis for given lot)
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	AF700

## Biological Properties

---

Appearance	Dark blue liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with AF700 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging
Recommended Dilutions	For flow cytometry applications, the suggested concentration is at 5 uL/million cells in 100 uL staining buffer. For the best performance of each application, the optimal concentration of this reagent needs to be carefully determined. <i>*The suggested working dilution is provided as a guide only. It is recommended that the users titrates the product for use in their tests using proper positive and negative controls.</i>

## Spectral Properties

---

Conjugate	AF700
-----------	-------

Excitation Wavelength 696 nm

Emission Wavelength 719 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 XFD700 (ex/em = 696/719 nm). XFD700 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 700 (Alexa Fluor® is the trademark of Thermo Fisher).