

iFluor™ 633 Anti-pig/ horse/ human/ non-human primates CD90 Antibody *5E10*Catalog number: 109000E0, 109000E1
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™ 633

Biological Properties

Appearance	Black liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 633 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 633
Excitation Wavelength	640 nm
Emission Wavelength	654 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™ 633 (ex/em = 640/654 nm). It is compatible with the 640 nm laser and 660/10 nm bandpass filter (for example, as in the BD FACSMelody™).