

**PE/iFluor™ 700 Anti-human CD34 Antibody
*4H11***Catalog number: 103401X0, 103401X1, 103401X2
Unit size: 25 tests, 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
Immunogen	CD34 (Gp105-120)
Clone	4H11
Conjugate	PE/iFluor™ 700

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PE/iFluor™ 700 under optimal conditions
Application	Flow Cytometry (FACS)

Spectral Properties

Conjugate	PE/iFluor™ 700
Excitation Wavelength	566 nm
Emission Wavelength	708 nm

Applications

4H11 is an anti-human monoclonal antibody that recognizes the CD34 antigen. CD34 (sometimes called Gp105-120) is a 105 - 120 kD glycoprotein that is located on the surface of cells like endothelial cells and stem cells. In many organisms, CD34 plays a role in the downregulation of tumor necrosis factor production, positively regulates granulocyte colony-stimulating factor production and is involved in the positive regulation of vasculogenesis. Also, it has been associated with vital biological processes such as endothelium development, especially glomerular endothelium development. From a research standpoint, it is of biological interest due to its association with critical

macromolecules/ligands like MadCAM-1, CRKL and L-Selectin. CD34 is a very popular antibody target, with over 55000 publications in the last decade. CD34 is vital to neuroscience, neuroinflammation and immunology research, typically serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to PE/iFluor™ 700 (ex/em = 566/708 nm). It is compatible with the 561 nm laser and 710/50 nm bandpass filter (for example, as in the BD LSRFortessa™ X-20).