

AF488 Anti-human CD34 Antibody *4H11*Catalog number: 10340150, 10340151
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	Human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Isotype	Mouse IgG1
Immunogen	CD34 (Gp105-120)
Clone	4H11
Conjugate	AF488

Biological Properties

Appearance	Red liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with AF488 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	AF488
Excitation Wavelength	499 nm
Emission Wavelength	520 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 AF488 (ex/em = 499/520 nm). It is compatible with the 488 nm laser and 530/30 nm bandpass filter (for example, as in the BD FACSCelesta™).