

FITC Anti-human CD135 Antibody *BV10A4*Catalog number: 113501G0, 113501G1
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	CD135 (FLT3, FLK2, STK-1)
Clone	BV10A4
Conjugate	FITC

Biological Properties

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

Spectral Properties

Conjugate	FITC
Excitation Wavelength	491 nm
Emission Wavelength	516 nm

Applications

BV10A4 is an anti-human monoclonal antibody that forms an immune complex with the CD135 antigen. CD135 (also known as FLT3, FLK2 or STK-1) is a 130 - 160 kD member of the immunoglobulin supergene family that is located on the surface of cells like stem cells. CD135 acts in important cellular pathways, in particular, the transmembrane receptor protein tyrosine kinase signaling pathway and cytokine-mediated signaling pathway. Also, in certain organisms, it positively regulates tyrosine phosphorylation of STAT protein, positively regulates phosphatidylinositol 3-kinase activity and is a promoter of phosphatidylinositol 3-kinase signaling. From a research standpoint, it is of biological interest due to its association with important macromolecules/ligands like FL (flt3 ligand). CD135 is a relatively rare antibody target, with fewer

than 900 publications in the last decade. Even still, CD135 has a variety of applications in immunology and cell biology research, frequently serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to FITC (ex/em = 491/516 nm). It is compatible with the 488 nm laser and 525/50 nm bandpass filter (for example, as in the Miltenyi Biotec MACSQuant X).