

iFluor™ 430 Anti-human CD135 Antibody
BV10A4Catalog number: 11350030, 11350031
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 | iFluor™ 430 |

Biological Properties

| | |
|-------------|--|
| Appearance | Yellow liquid |
| Preparation | Antibody purified by affinity chromatography and then conjugated with iFluor™ 430 under optimal conditions |
| Application | Flow Cytometry (FACS), Fluorescence Imaging |

Spectral Properties

| | |
|-----------------------|-------------|
| Conjugate | iFluor™ 430 |
| Excitation Wavelength | 433 nm |
| Emission Wavelength | 498 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 iFluor™ 430 (ex/em = 433/498 nm). It is compatible with the 445 nm laser and 510/80 nm bandpass filter (for example, as in the BD FACS Aria™ III).