

**mFluor™ Violet 450 Anti-human CD151  
Antibody \*50-6\***Catalog number: 115100Z0, 115100Z1  
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 kappa            |
| Immunogen          | CD151 (PETA-3, SFA-1, GP27) |
| Clone              | 50-6                        |
| Conjugate          | mFluor™ Violet 450          |

**Biological Properties**

|             |                                                                                                                   |
|-------------|-------------------------------------------------------------------------------------------------------------------|
| Appearance  | Light yellow liquid                                                                                               |
| Preparation | Antibody purified by affinity chromatography and then conjugated with mFluor™ Violet 450 under optimal conditions |
| Application | Flow Cytometry (FACS), Fluorescence Imaging                                                                       |

**Spectral Properties**

|                       |                    |
|-----------------------|--------------------|
| Conjugate             | mFluor™ Violet 450 |
| Excitation Wavelength | 406 nm             |
| Emission Wavelength   | 445 nm             |

**Applications**

50-6 is an anti-human monoclonal antibody that is specific for the CD151 antigen. CD151 (alternatively called PETA-3, SFA-1 or GP27) is a 29 kD member of the tetraspanin family that is located on the surface of cells such as stem cells, epithelial cells, endothelial cells and platelets. CD151 is associated with a variety of biologically interesting macromolecules/ligands, namely,  $\alpha 6$  and integrins  $\alpha 3$ . CD151 is a fairly uncommon

antibody target, with a little more than 1600 publications in the last decade. Even still, CD151 is vital to stem cells 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 mFluor™ Violet 450 (ex/em = 406/445 nm). It is compatible with the 405 nm laser and 450/50 nm bandpass filter (for example, as in the BD FACSaria™ Fusion).