

**mFluor™ Violet 510 Anti-human CD49  
Antibody \*ASC-1\***Catalog number: 10492110, 10492111  
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          | CD49c (ITGA3)      |
| Clone              | ASC-1              |
| Conjugate          | mFluor™ Violet 510 |

**Biological Properties**

|             |                                                                                                                   |
|-------------|-------------------------------------------------------------------------------------------------------------------|
| Appearance  | Yellow liquid                                                                                                     |
| Preparation | Antibody purified by affinity chromatography and then conjugated with mFluor™ Violet 510 under optimal conditions |
| Application | Flow Cytometry (FACS), Fluorescence Imaging                                                                       |

**Spectral Properties**

|                       |                    |
|-----------------------|--------------------|
| Conjugate             | mFluor™ Violet 510 |
| Excitation Wavelength | 412 nm             |
| Emission Wavelength   | 505 nm             |

**Applications**

ASC-1 is an anti-human monoclonal antibody that targets the CD49c antigen. CD49c (sometimes called ITGA3) is a 150 kD member of the integrin family that is expressed on the surface of cells like B cells, T cells and epithelial cells. CD49 is associated with a variety of biologically interesting macromolecules/ligands, for instance, laminin, CD9 and collagen. CD49 is a relatively rare antibody target, with fewer than 600

publications in the last decade. Even still, CD49c has a variety of applications in neuroscience and immunology research, commonly 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 510 (ex/em = 412/505 nm). It is compatible with the 405 nm laser and 528/45 nm bandpass filter (for example, as in the BD FACSMelody™).