

**iFluor™ 840 Anti-human CD49 Antibody**  
**\*ASC-1\***Catalog number: 104920Q0, 104920Q1  
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          | iFluor™ 840      |

**Biological Properties**

|             |                                                                                                            |
|-------------|------------------------------------------------------------------------------------------------------------|
| Appearance  | Dark brown liquid                                                                                          |
| Preparation | Antibody purified by affinity chromatography and then conjugated with iFluor™ 840 under optimal conditions |
| Application | Flow Cytometry (FACS), Fluorescence Imaging                                                                |

**Spectral Properties**

|                       |             |
|-----------------------|-------------|
| Conjugate             | iFluor™ 840 |
| Excitation Wavelength | 836 nm      |
| Emission Wavelength   | 879 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 iFluor™ 840 (ex/em = 836/879 nm).