

**APC/iFluor™ 700 Anti-human CD49
Antibody *ASC-1***Catalog number: 104921F0, 104921F1, 104921F2
Unit size: 25 tests, 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	APC/iFluor™ 700

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with APC/iFluor™ 700 under optimal conditions
Application	Flow Cytometry (FACS)

Spectral Properties

Conjugate	APC/iFluor™ 700
Excitation Wavelength	685 nm
Emission Wavelength	710 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 APC/iFluor™ 700 (ex/em = 685/710 nm).