

**APC/Cy7 Anti-mouse CD49 Antibody  
\*5H10-27 (MFR5)\***Catalog number: 104931D0, 104931D1, 104931D2  
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 | Mouse                                                     |
| Class              | Primary                                                   |
| Clonality          | Monoclonal                                                |
| Host               | Rat                                                       |
| Isotype            | Rat IgG2a kappa                                           |
| Immunogen          | CD49e (VLA-5 $\alpha$ , Integrin $\alpha$ 5 chain, ITGA5) |
| Clone              | 5H10-27 (MFR5)                                            |
| Conjugate          | APC/Cy7                                                   |

**Biological Properties**

|             |                                                                                                        |
|-------------|--------------------------------------------------------------------------------------------------------|
| Preparation | Antibody purified by affinity chromatography and then conjugated with APC/Cy7 under optimal conditions |
| Application | Flow Cytometry (FACS)                                                                                  |

**Spectral Properties**

|                       |         |
|-----------------------|---------|
| Conjugate             | APC/Cy7 |
| Excitation Wavelength | 754 nm  |
| Emission Wavelength   | 779 nm  |

**Applications**

5H10-27 (MFR5) is an anti-mouse monoclonal antibody that forms an immune complex with the CD49e antigen. CD49e (also known as ITGA5, Integrin  $\alpha$ 5 chain or VLA-5 $\alpha$ ) is a 135 kD member of the Integrin family that is found on the surface of cells like stem cells and T cells. CD49 is associated with a variety of biologically interesting macromolecules/ligands, for example, fibronectin. CD49 is a relatively rare antibody target, with fewer than 600 publications in the last decade. Even still, CD49e is commonly used in flow cytometry applications as a phenotypic marker for differentiation of cell types, especially in the study of cell biology. This antibody was purified through affinity chromatography and conjugated to APC/Cy7 (ex/em = 754/779 nm).