

**PerCP/Cy5.5 Anti-mouse/human CD49d  
Antibody \*PS/2\***Catalog number: 104901U0, 104901U1, 104901U2  
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, human                                        |
| Class              | Primary                                             |
| Clonality          | Monoclonal                                          |
| Host               | Rat                                                 |
| Isotype            | Rat igg2b, $\kappa$                                 |
| Immunogen          | CD49d (VLA-4 $\alpha$ , ITGA4, Integrin $\alpha$ 4) |
| Clone              | PS/2                                                |
| Conjugate          | PerCP/Cy5.5                                         |

**Biological Properties**

|             |                                                                                                            |
|-------------|------------------------------------------------------------------------------------------------------------|
| Preparation | Antibody purified by affinity chromatography and then conjugated with PerCP/Cy5.5 under optimal conditions |
| Application | Flow Cytometry (FACS)                                                                                      |

**Spectral Properties**

|                       |             |
|-----------------------|-------------|
| Conjugate             | PerCP/Cy5.5 |
| Excitation Wavelength | 489 nm      |
| Emission Wavelength   | 679 nm      |

**Applications**

PS/2 is an anti-mouse/human monoclonal antibody that forms an immune complex with the CD49d antigen. CD49d (sometimes called  $\alpha$ 4 integrin, Integrin  $\alpha$ 4 chain, VLA-4  $\alpha$  chain or Integrin alpha-4) is a 150 kD glycoprotein that is found on the surface of cells such as NK cells, dendritic cells and macrophages. CD49d is associated with a variety of biologically interesting macromolecules/ligands, for instance, Paxillin and CD106. CD49d is a fairly uncommon antibody target, with a little more than 3400 publications in the last decade. Even still, CD49d is vital to cell biology and immunology 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 PerCP/Cy5.5 (ex/em = 489/679 nm). It is compatible with the 488 nm laser and 695/40 nm bandpass filter (for example, as in the BD FACS Aria™ III).