

**PE/Cy5 Anti-human CD328 Antibody**  
**\*6-434\***Catalog number: 132801N0, 132801N1, 132801N2  
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             |
| Immunogen          | CD328 (Siglec7, AIRM1) |
| Clone              | 6-434                  |
| Conjugate          | PE/Cy5                 |

**Biological Properties**

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

**Spectral Properties**

|                       |        |
|-----------------------|--------|
| Conjugate             | PE/Cy5 |
| Excitation Wavelength | 565 nm |
| Emission Wavelength   | 666 nm |

**Applications**

6-434 is an anti-human monoclonal antibody that forms an immune complex with the CD328 antigen. CD328 (alternatively called Siglec7 or AIRM1) is a 75 kD transmembrane protein that is found on the surface of cells like macrophages and T cells. CD328 is associated with a variety of biologically interesting macromolecules/ligands, in particular, silylated glycans. CD328 is a relatively rare antibody target, with fewer than 20 publications in the last decade. Even still, CD328 has a variety of applications in immunology research, often serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to PE/Cy5 (ex/em = 565/666 nm). It is compatible with the 561 nm laser and 660/20 nm bandpass filter (for example, as in the Agilent Technologies

NovoCyte).