

**PacOrange Anti-human CD235 Antibody**  
**\*HI264\***Catalog number: 123501K0, 123501K1  
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 IgG2a                                   |
| Immunogen          | CD235a (GYPA, Sialoglycoprotein alpha, PAS-2) |
| Clone              | HI264                                         |
| Conjugate          | PacOrange                                     |

**Biological Properties**

|             |                                                                                                          |
|-------------|----------------------------------------------------------------------------------------------------------|
| Preparation | Antibody purified by affinity chromatography and then conjugated with PacOrange under optimal conditions |
| Application | Flow Cytometry (FACS), Fluorescence Imaging                                                              |

**Spectral Properties**

|                       |           |
|-----------------------|-----------|
| Conjugate             | PacOrange |
| Excitation Wavelength | 400 nm    |
| Emission Wavelength   | 551 nm    |

**Applications**

HI264 is an anti-human monoclonal antibody that forms an immune complex with the CD235a antigen. CD235a (sometimes called MN sialoglycoprotein or MNS blood group) is a 10 kD single-pass type I membrane protein that is expressed on the surface of cells like erythrocytes. CD235 is associated with a variety of biologically interesting macromolecules/ligands, namely, Plasmodium falciparum erythrocyte binding antigen EBA-175, CD170 and Influenza virus. CD235 is a relatively rare antibody target, with fewer than 6000 publications in the last decade. Even still, CD235a is essential for cell adhesion and 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 PacOrange (ex/em = 400/551 nm). It is compatible with the 405 nm laser and 525/50 nm bandpass filter (for example, as in the Miltenyi Biotec MACSQuant X).