

**PacBlue Anti-human CD172a Antibody**  
**\*15-414\***Catalog number: 11720110, 11720111  
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          | CD172a (SHPS-1, SIRP $\alpha$ , BIT, MFR, MYD-1, P84) |
| Clone              | 15-414                                                |
| Conjugate          | PacBlue                                               |

**Biological Properties**

---

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

**Spectral Properties**

---

|                       |         |
|-----------------------|---------|
| Conjugate             | PacBlue |
| Excitation Wavelength | 404 nm  |
| Emission Wavelength   | 455 nm  |

**Applications**

---

The 15-414 monoclonal antibody binds with human CD172a, a 85 - 90 kD glycoprotein often found on the surface of dendritic cells and neurons. CD172a is associated with a variety of biologically interesting macromolecules/ligands, in particular, CD47 and PTPN11. CD172a is a relatively rare antibody target, with fewer than 700 publications in the last decade. Even still, CD172a has a variety of applications in immunology and cell

biology 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 PacBlue (ex/em = 404/455 nm). It is compatible with the 405 nm laser and 450/50 nm bandpass filter (for example, as in the Miltenyi Biotec MACSQuant Analyzer 16).