

**PE/Texas Red® Anti-human CD111 Antibody**  
**\*R1.302\***Catalog number: 111101R0, 111101R1, 111101R2  
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 kappa                    |
| Immunogen          | CD111 (Nectin-1, HVEC, PVRL1, PRR1) |
| Clone              | R1.302                              |
| Conjugate          | PE/Texas Red®                       |

**Biological Properties**

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

**Spectral Properties**

|                       |               |
|-----------------------|---------------|
| Conjugate             | PE/Texas Red® |
| Excitation Wavelength | 567 nm        |
| Emission Wavelength   | 615 nm        |

**Applications**

R1.302 is an anti-human monoclonal antibody that forms an immune complex with the CD111 antigen. CD111 (also known as PVRL1 or Nectin-1) is a 75 kD single-pass type I membrane protein that is found on the surface of cells like erythrocytes, endothelial cells, epithelial cells, stem cells and macrophages. CD111 has been thought to be involved with critical biological processes like cell-cell adhesion, particularly homophilic cell adhesion via plasma membrane adhesion molecules, and is associated with a variety of biologically interesting macromolecules/ligands, for instance, nectin3 and afadin gd. CD111 is a relatively rare antibody target, with fewer than 90 publications in the

last decade. Even still, CD111 is often used in flow cytometry applications as a phenotypic marker for differentiation of cell types, specifically in the study of neuroscience. This antibody was purified through affinity chromatography and conjugated to PE/Texas Red® (ex/em = 567/615 nm). It is compatible with the 561 nm laser and 615/24 nm bandpass filter (for example, as in the Bio-Rad ZES Cell Analyzer).