

**PE/Texas Red® Anti-human CD102 Antibody**  
**\*CBR-IC2/2\***Catalog number: 110201R0, 110201R1, 110201R2  
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 IgG2a   |
| Immunogen          | CD102 (ICAM2) |
| Clone              | CBR-IC2/2     |
| 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**

CBR-IC2/2 is an anti-human monoclonal antibody that targets the CD102 antigen. CD102 (alternatively called ICAM2) is a 55 kD transmembrane glycoprotein that is expressed on the surface of cells like T cells and macrophages. CD102 is associated with a variety of biologically interesting macromolecules/ligands, for instance, LFA-1, integrin  $\alpha\beta 2$  and CD11b/CD18. CD102 is a relatively rare antibody target, with fewer than 200 publications in the last decade. Even still, CD102 is often used in flow cytometry applications as a phenotypic marker for differentiation of cell types, specifically in the study of immunology. 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 615/20 nm bandpass filter (for example, as in the Agilent Technologies NovoCyte).