

**APC/Cy7 Anti-human CD19 Antibody \*4G7\***Catalog number: 101931D0, 101931D1, 101931D2  
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          | CD19 (B4)  |
| Clone              | 4G7        |
| Conjugate          | APC/Cy7    |

**Biological Properties**

|             |                                                                                                        |
|-------------|--------------------------------------------------------------------------------------------------------|
| Preparation | Antibody purified by affinity chromatography and then conjugated with APC/Cy7 under optimal conditions |
| Application | Flow Cytometry (FACS)                                                                                  |

**Spectral Properties**

|                       |         |
|-----------------------|---------|
| Conjugate             | APC/Cy7 |
| Excitation Wavelength | 754 nm  |
| Emission Wavelength   | 779 nm  |

**Applications**

The 4G7 monoclonal antibody binds with human CD19, a 95 kD transmembrane protein commonly expressed on the surface of B cells and lymphocytes. CD19 plays a role in essential cellular pathways, for example, the antigen receptor-mediated signaling pathway and B cell receptor signaling pathway. Moreover, in some organisms, it acts to positively regulate phosphatidylinositol 3-kinase activity, is an enhancer of protein kinase B signaling and is involved in the positive regulation of release of sequestered calcium ion into cytosol. From a research standpoint, it is of biological interest due to its association with important macromolecules/ligands such as Fyn and PI3-kinase. CD19 is a very popular antibody target, with over 30000 publications in the last decade. CD19 is essential for immunology research, commonly serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to

APC/Cy7 (ex/em = 754/779 nm).