

**FITC Anti-human/monkey CD20 Antibody**  
**\*2H7\***Catalog number: 102021G0, 102021G1  
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, monkey   |
| Class              | Primary         |
| Clonality          | Monoclonal      |
| Host               | Mouse           |
| Isotype            | Mouse igg2b, κ  |
| Immunogen          | CD20 (Bp35, B1) |
| Clone              | 2H7             |
| Conjugate          | FITC            |

**Biological Properties**

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

**Spectral Properties**

|                       |        |
|-----------------------|--------|
| Conjugate             | FITC   |
| Excitation Wavelength | 491 nm |
| Emission Wavelength   | 516 nm |

**Applications**

The 2H7 monoclonal antibody binds with human/monkey CD20, a 33 - 37 kD transmembrane glycoprotein commonly expressed on the surface of T cells and B cells. CD20 is a component of essential cellular pathways, in particular, the B cell receptor signaling pathway. From a research standpoint, it is of biological interest due to its association with vital macromolecules/ligands such as Lyn. CD20 is a moderately popular antibody target, with over 20000 publications in the last decade. CD20 is vital to immunology 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 FITC (ex/em = 491/516 nm). It is compatible with the 488 nm laser and 530/30 nm bandpass filter (for example, as in the BD FACS Aria™ II).