

**mFluor™ Green 620 Anti-human/monkey  
CD20 Antibody \*2H7\***Catalog number: 102020U0, 102020U1  
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          | mFluor™ Green 620 |

**Biological Properties**

|             |                                                                                                                  |
|-------------|------------------------------------------------------------------------------------------------------------------|
| Appearance  | Purple liquid                                                                                                    |
| Preparation | Antibody purified by affinity chromatography and then conjugated with mFluor™ Green 620 under optimal conditions |
| Application | Flow Cytometry (FACS), Fluorescence Imaging                                                                      |

**Spectral Properties**

|                       |                   |
|-----------------------|-------------------|
| Conjugate             | mFluor™ Green 620 |
| Excitation Wavelength | 525 nm            |
| Emission Wavelength   | 623 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 mFluor™ Green 620 (ex/em = 525/623 nm). It is compatible with the 532 nm laser and 609/30 nm bandpass filter (for example, as in the Luminex Guava easyCyte).