

**mFluor™ Green 620 Anti-human CD20
Antibody *HI20a***Catalog number: 102010U0, 102010U1
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 |
| Class | Primary |
| Clonality | Monoclonal |
| Host | Mouse |
| Isotype | Mouse IgG2a |
| Immunogen | CD20 (Bp35, B1) |
| Clone | HI20a |
| 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

HI20a is an anti-human monoclonal antibody that recognizes the CD20 antigen. CD20 (alternatively called MS4A1, Bp35 or Leukocyte surface antigen Leu-16) is a 33 - 37 kD transmembrane protein that is expressed on the surface of cells such as B cells. CD20 plays a role in important cellular pathways, namely, the B cell receptor signaling pathway. From a research standpoint, it is of biological interest due to its association

with key macromolecules/ligands such as Lyn. CD20 is a moderately popular antibody target, with over 20000 publications in the last decade. CD20 has a variety of applications in costimulatory molecules and 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 mFluor™ Green 620 (ex/em = 525/623 nm). It is compatible with the 532 nm laser and 620/15 nm bandpass filter (for example, as in the Thermo Fisher Attune NxT).