

**PE/Texas Red® Anti-human CD20 Antibody
*HI47***Catalog number: 102001S0, 102001S1, 102001S2
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 IgG3 |
| Immunogen | CD20 (Bp35, B1) |
| Clone | HI47 |
| 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

HI47 is an anti-human monoclonal antibody that recognizes the CD20 antigen. CD20 (alternatively called Leukocyte surface antigen Leu-16) is a 33 - 37 kD transmembrane protein that is located on the surface of cells such as B cells and T cells. CD20 is a member of critical cellular pathways, in particular, the B cell receptor signaling pathway. From a research standpoint, it is of biological interest due to its association with essential macromolecules/ligands like lyn. CD20 is a very popular antibody target, with over 28000 publications in the last decade. CD20 has been widely used in immunology and costimulatory molecules research, often serving as a phenotypic marker for differentiating cell types in

flow cytometric applications. 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 615/24 nm bandpass filter (for example, as in the Bio-Rad ZE5 Cell Analyzer).