

**PE/Cy5 Anti-human CD52 Antibody \*HI186\***Catalog number: 105201M0, 105201M1, 105201M2  
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 IgG2b           |
| Immunogen          | CD52 (CAMPATH-1, HE5) |
| Clone              | HI186                 |
| Conjugate          | PE/Cy5                |

**Biological Properties**

---

|             |   |
|-------------|---|
| Preparation | Antibody purified by affinity chromatography and then conjugated with PE/Cy5 under optimal conditions |
| Application | Flow Cytometry (FACS)   |

**Spectral Properties**

---

|                       |        |
|-----------------------|--------|
| Conjugate             | PE/Cy5 |
| Excitation Wavelength | 565 nm |
| Emission Wavelength   | 666 nm |

**Applications**

---

HI186 is an anti-human monoclonal antibody that is specific for the CD52 antigen. CD52 (also known as HE5) is a 7 kD transmembrane glycoprotein that is located on the surface of cells such as epithelial cells, NK cells and T cells. In some organisms, CD52 is an enhancer of cytosolic calcium ion concentration, and is associated with a variety of biologically interesting macromolecules/ligands, in particular, . CD52 is a fairly uncommon antibody target, with a little more than 2600 publications in the last decade. Even still, CD52 is vital to 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/Cy5 (ex/em = 565/666 nm). It is compatible with the 561 nm laser and 670/30 nm bandpass filter (for example, as in the BD LSRFortessa™ X-20).