

**PerCP Anti-human CD58 Antibody \*HI58a\***Catalog number: 105801T0, 105801T1, 105801T2  
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 IgG1   |
| Immunogen          | CD58 (LFA-3) |
| Clone              | HI58a        |
| Conjugate          | PerCP        |

**Biological Properties**

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

**Spectral Properties**

|                       |        |
|-----------------------|--------|
| Conjugate             | PerCP  |
| Excitation Wavelength | 477 nm |
| Emission Wavelength   | 678 nm |

**Applications**

HI58a is an anti-human monoclonal antibody that targets the CD58 antigen. CD58 (sometimes referred to as Lymphocyte function-associated antigen 3, LFA3 or LFA-3) is a 45 - 70 kD transmembrane protein that is located on the surface of cells such as macrophages, NK cells, dendritic cells, B cells and granulocytes. CD58 has been associated with key biological processes such as cell-cell adhesion, especially heterotypic cell-cell adhesion. Also, in certain organisms, it is involved in the positive regulation of interleukin-8 secretion. From a research standpoint, it is of biological interest due to its association with important macromolecules/ligands like CD2 and LFA-2. CD58 is a relatively rare antibody target, with less than 1000 publications in the last decade. Even still, CD58 is commonly used in flow cytometry applications as a phenotypic marker for differentiation of cell types, especially in the study of cell biology. This antibody was purified through affinity chromatography and conjugated to

PerCP (ex/em = 477/678 nm). It is compatible with the 488 nm laser and 700/54 nm bandpass filter (for example, as in the BD FACSVerse™).