

**iFluor™ 568 Anti-human CD62L Antibody**  
**\*HI62L\***Catalog number: 106210B0, 106210B1  
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          | CD62L (L-selectin, LECAM-1, LAM-1, Leu-8, TQ-1) |
| Clone              | HI62L   |
| Conjugate          | iFluor™ 568                                     |

**Biological Properties**

|             |  |
|-------------|--|
| Appearance  | Purple liquid  |
| Preparation | Antibody purified by affinity chromatography and then conjugated with iFluor™ 568 under optimal conditions |
| Application | Flow Cytometry (FACS), Fluorescence Imaging  |

**Spectral Properties**

|                       |             |
|-----------------------|-------------|
| Conjugate             | iFluor™ 568 |
| Excitation Wavelength | 568 nm      |
| Emission Wavelength   | 587 nm      |

**Applications**

The HI62L monoclonal antibody binds to human CD62L, a 74 - 95 kD single-pass type I membrane protein typically located on the surface of T cells, monocytes, neutrophils and thymocytes. CD62L is associated with a variety of biologically interesting macromolecules/ligands, for example, MADCAM-1. CD62L is a fairly uncommon antibody target, with a little more than 10000 publications in the last decade. Even still, CD62L

has been widely used in innate immunity 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 iFluor™ 568 (ex/em = 568/587 nm). It is compatible with the 561 nm laser and 577/35 nm bandpass filter (for example, as in the Luminex Amnis ImageStream).