

**iFluor™ 700 Anti-human CD62I Antibody**  
**\*HI62L\***Catalog number: 106210J0, 106210J1  
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	CD62I (L-selectin, LECAM-1, LAM-1, Leu-8, TQ-1)
Clone	HI62L
Conjugate	iFluor™ 700

**Biological Properties**

---

Appearance	Blue liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 700 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

**Spectral Properties**

---

Conjugate	iFluor™ 700
Excitation Wavelength	690 nm
Emission Wavelength	713 nm

**Applications**

---

The HI62L monoclonal antibody binds to human CD62I, a 74 - 95 kD single-pass type I membrane protein typically located on the surface of T cells, monocytes, neutrophils and thymocytes. CD62I is associated with a variety of biologically interesting macromolecules/ligands, for example, MADCAM-1. CD62I is a fairly uncommon antibody target, with a little more than 10000 publications in the last decade. Even still, CD62I

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™ 700 (ex/em = 690/713 nm).