

**mFluor™ Red 700 Anti-human CD62  
Antibody \*HI62E\***Catalog number: 106200V0, 106200V1  
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	CD62e (E-selectin, ELAM-1)
Clone	HI62E
Conjugate	mFluor™ Red 700

**Biological Properties**

---

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

**Spectral Properties**

---

Conjugate	mFluor™ Red 700
Excitation Wavelength	680 nm
Emission Wavelength	695 nm

**Applications**

---

The HI62E monoclonal antibody reacts with human CD62e, a 115 kD single-pass type I membrane protein frequently expressed on the surface of endothelial cells and  $\text{tnf } \alpha$ s. CD62 has been associated with vital biological processes like inflammatory response, particularly leukocyte migration involved in inflammatory response. Also, in many organisms, it is an enhancer of receptor internalization. From a research standpoint,

it is of biological interest due to its association with key macromolecules/ligands such as a and Sialyl Lewis x. CD62 is a relatively rare antibody target, with fewer than FAKE103000FAKE publications in the last decade. Even still, CD62e is often used in flow cytometry applications as a phenotypic marker for differentiation of cell types, specifically in the study of cell biology, neuroscience cell markers and immunology. This antibody was purified through affinity chromatography and conjugated to mFluor™ Red 700 (ex/em = 680/695 nm). It is compatible with the 642 nm laser and 702/87 nm bandpass filter (for example, as in the Luminex Amnis CellStream).