

PE/iFluor™ 750 Anti-human CD62L Antibody
HI62LCatalog number: 106211R0, 106211R1, 106211R2
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 IgG2a
Immunogen	CD62L (L-selectin, LECAM-1, LAM-1, Leu-8, TQ-1)
Clone	HI62L
Conjugate	PE/iFluor™ 750

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PE/iFluor™ 750 under optimal conditions
Application	Flow Cytometry (FACS)

Spectral Properties

Conjugate	PE/iFluor™ 750
Excitation Wavelength	566 nm
Emission Wavelength	778 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 PE/iFluor™ 750 (ex/em = 566/778 nm). It is

compatible with the 561 nm laser and 780/60 nm bandpass filter (for example, as in the BD FACSCelesta™).