

PE/XFD610 Anti-human CD62I Antibody *HI62L*

Catalog Number: 106211O0,
106211O1, 106211O2
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	Lot specific (please consult certificate of analysis for given lot)
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	PE/AF610

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PE/AF610 under optimal conditions
Application	Flow Cytometry (FACS)
Recommended Dilutions	For flow cytometry applications, the suggested concentration is at 5 uL/million cells in 100 uL staining buffer. For the best performance of each application, the optimal concentration of this reagent needs to be carefully determined.
	<i>*The suggested working dilution is provided as a guide only. It is recommended that the users titrates the product for use in their tests using proper positive and negative controls.</i>

Spectral Properties

Conjugate	PE/AF610
-----------	----------

Excitation Wavelength 565 nm

Emission Wavelength 627 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/XFD610 (ex/em = 567/627 nm). XFD610 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 610 (Alexa Fluor® is the trademark of Thermo Fisher). It is compatible with the 561 nm laser and 610/30 nm bandpass filter (for example, as in the Luminex Amnis FlowSight).