

PacOrange Anti-human CD62 Antibody
HI62ECatalog number: 106201K0, 106201K1
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	PacOrange

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PacOrange under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	PacOrange
Excitation Wavelength	400 nm
Emission Wavelength	551 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 103,000 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 PacOrange (ex/em = 400/551 nm). It is compatible with the 405 nm laser and 525/50 nm bandpass filter (for example, as in the BD LSRFortessa™ Cell Analyzer).