

**APC/iFluor™ 700 Anti-human CD62
Antibody *HI62E***Catalog number: 106201F0, 106201F1, 106201F2
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	CD62e (E-selectin, ELAM-1)
Clone	HI62E
Conjugate	APC/iFluor™ 700

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with APC/iFluor™ 700 under optimal conditions
Application	Flow Cytometry (FACS)

Spectral Properties

Conjugate	APC/iFluor™ 700
Excitation Wavelength	685 nm
Emission Wavelength	710 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 APC/iFluor™ 700 (ex/em = 685/710 nm).