

**PacOrange Anti-mouse CD106 Antibody
*429 (MVCAM.A)***Catalog number: 110601J0, 110601J1
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	Mouse
Class	Primary
Clonality	Monoclonal
Host	Rat
Isotype	Rat IgG2a kappa
Immunogen	CD106 (VCAM1, INCAM-110)
Clone	429 (MVCAM.A)
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 429 (MVCAM.A) monoclonal antibody binds with mouse CD106, a 110 kD transmembrane protein frequently found on the surface of dendritic cells and endothelial progenitors. CD106 is associated with a variety of biologically interesting macromolecules/ligands, for instance, integrin $\alpha 4\beta 1$ and VLA-4. CD106 is a fairly uncommon antibody target, with a little more than 2300 publications in the last decade. Even still, CD106 has a variety of applications in stem cells and cell adhesion research, often serving as a phenotypic marker for differentiating cell types in flow cytometric applications. 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 Bio-Rad ZE5 Cell Analyzer).