

**iFluor™ 647 Anti-mouse CD106 Antibody
*429 (MVCAM.A)***Catalog number: 110600F0, 110600F1
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	iFluor™ 647

Biological Properties

Appearance	Blue liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 647 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 647
Excitation Wavelength	656 nm
Emission Wavelength	670 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 $\alpha\beta 1$ 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 iFluor™ 647 (ex/em = 656/670 nm). It is compatible with the 640 nm laser and 660/20 nm bandpass filter (for example, as in the BD FACS Aria™ II).