

**iFluor™ 710 Anti-mouse CD106 Antibody  
\*429 (MVCAM.A)\***Catalog number: 110600K0, 110600K1  
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™ 710              |

**Biological Properties**

|             |  |
|-------------|--|
| Appearance  | Blue liquid  |
| Preparation | Antibody purified by affinity chromatography and then conjugated with iFluor™ 710 under optimal conditions |
| Application | Flow Cytometry (FACS), Fluorescence Imaging  |

**Spectral Properties**

|                       |             |
|-----------------------|-------------|
| Conjugate             | iFluor™ 710 |
| Excitation Wavelength | 717 nm      |
| Emission Wavelength   | 739 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™ 710 (ex/em = 717/739 nm).