

**Purified Anti-mouse CD106 Antibody *429
(MVCAM.A)***

Catalog number: 11060000

Unit size: 100 ug

Product Details

Storage Conditions	2-8°C with minimized light exposure. Do not freeze.
Expiration Date	12 months upon receiving
Concentration	0.5 mg/mL
Formulation	Phosphate-buffered saline (PBS, pH 7.2), 0.09% sodium azide

Antibody Properties

Species Reactivity	Mouse
Class	Primary
Clonality	Monoclonal
Host	Rat
Isotype	Rat IgG2a kappa
Immunogen	CD106 (VCAM1, INCAM-110)
Clone	429 (MVCAM.A)

Biological Properties

Appearance	Liquid
Preparation	Antibody purified by affinity chromatography
Application	Flow Cytometry (FACS), ELISA, HC, Western Blot

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.