

**iFluor™ 860 Anti-mouse/human CD49d
Antibody *PS/2***Catalog number: 104900R0, 104900R1
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, human
Class	Primary
Clonality	Monoclonal
Host	Rat
Isotype	Rat igg2b, κ
Immunogen	CD49d (VLA-4 α , ITGA4, Integrin α 4)
Clone	PS/2
Conjugate	iFluor™ 860

Biological Properties

Appearance	Green liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 860 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 860
Excitation Wavelength	853 nm
Emission Wavelength	878 nm

Applications

PS/2 is an anti-mouse/human monoclonal antibody that forms an immune complex with the CD49d antigen. CD49d (sometimes called α 4 integrin, Integrin α 4 chain, VLA-4 α chain or Integrin alpha-4) is a 150 kD glycoprotein that is found on the surface of cells such as NK cells, dendritic cells and macrophages. CD49d is associated with a variety of biologically interesting macromolecules/ligands, for instance, Paxillin and

CD106. CD49d is a fairly uncommon antibody target, with a little more than 3400 publications in the last decade. Even still, CD49d is vital to cell biology and immunology research, frequently serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to iFluor™ 860 (ex/em = 853/878 nm).