

PE/XFD700 Anti-mouse/human CD49d Antibody *PS/2*

Catalog Number: 104901P0,
104901P1, 104901P2
Unit Size: 25 tests, 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	Lot specific (please consult certificate of analysis for given lot)
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	PE/AF700

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PE/AF700 under optimal conditions
Application	Flow Cytometry (FACS)
Recommended Dilutions	For flow cytometry applications, the suggested concentration is at 5 μ L/million cells in 100 μ L staining buffer. For the best performance of each application, the optimal concentration of this reagent needs to be carefully determined.
	<i>*The suggested working dilution is provided as a guide only. It is recommended that the users titrates the product for use in their tests using proper positive and negative controls.</i>

Spectral Properties

Conjugate	PE/AF700
-----------	----------

Excitation Wavelength 565 nm

Emission Wavelength 721 nm

Applications

PS/2 is an anti-mouse/human monoclonal antibody that forms an immune complex with the CD49d antigen. CD49d (sometimes called $\alpha 4$ integrin, Integrin $\alpha 4$ 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 PE/XFD700 (ex/em = 566/721 nm). XFD700 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 700 (Alexa Fluor® is the trademark of Thermo Fisher). It is compatible with the 561 nm laser and 702/87 nm bandpass filter (for example, as in the Luminex Amnis CellStream).