

XFD680 Anti-mouse/human CD49d Antibody *PS/2*

Catalog Number: 10490190, 10490191

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	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	AF680

Biological Properties

Appearance	Blue liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with AF680 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging
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	AF680
Excitation Wavelength	681 nm
Emission Wavelength	704 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 XFD680 (ex/em = 681/704 nm). XFD680 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 680 (Alexa Fluor® is the trademark of Thermo Fisher). It is compatible with the 642 nm laser and 702/87 nm bandpass filter (for example, as in the Luminex Amnis CellStream).