

APC/XFD750 Anti-human CD9 Antibody
***HI9a, XFD750 Same Structure to Alexa**
Fluor™ 750*Catalog number: 100901E0, 100901E1, 100901E2
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	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	Human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Isotype	Mouse IgG1
Immunogen	CD9 (MRP-1, P24)
Clone	HI9a
Conjugate	APC/AF750

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with APC/AF750 under optimal conditions
Application	Flow Cytometry (FACS)

Spectral Properties

Conjugate	APC/AF750
Excitation Wavelength	756 nm
Emission Wavelength	785 nm

Applications

HI9a is an anti-human monoclonal antibody that recognizes the CD9 antigen. CD9 (alternatively called MIC3, TSPAN29, Tetraspanin or 5H9 antigen) is a 24 kD member of the Tetraspan family that is expressed on the surface of cells like macrophages, granulocytes, epithelial cells, platelets and endothelial cells. In many organisms, CD9 is involved in the negative regulation of cellular component movement, represses cell population proliferation and negatively regulates platelet aggregation. From a research standpoint, it is of biological interest due to its

association with critical macromolecules/ligands such as Integrin β 1. CD9 is a fairly uncommon antibody target, with a little more than 6100 publications in the last decade. Even still, CD9 is frequently used in flow cytometry applications as a phenotypic marker for differentiation of cell types, particularly in the study of immunology and stem cells. This antibody was purified through affinity chromatography and conjugated to APC/XFD750 (ex/em = 756/785 nm). XFD750 is manufactured by AAT Bioquest, and it has the same chemical structure of Alexa Fluor® 750 (Alexa Fluor® is the trademark of ThermoFisher).