

APC Anti-human CD144 Antibody *55-7H1*Catalog number: 114401B0, 114401B1, 114401B2
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 kappa
Immunogen	CD144 (VE cadherin, cadherin-5)
Clone	55-7H1
Conjugate	APC

Biological Properties

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

Spectral Properties

Conjugate	APC
Excitation Wavelength	651 nm
Emission Wavelength	660 nm

Applications

55-7H1 is an anti-human monoclonal antibody that targets the CD144 antigen. CD144 (also known as VE cadherin or cadherin-5) is a member of the cadherin family that is expressed on the surface of cells like endothelial cells. CD144 is associated with a variety of biologically interesting macromolecules/ligands, in particular, plakoglobin. CD144 is a fairly uncommon antibody target, with a little more than 1500 publications in the last decade. Even still, CD144 is essential for cell motility/cytoskeleton/structure, cell biology and angiogenesis 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 APC (ex/em = 651/660 nm). It is compatible with the 642 nm laser and 664/20 nm bandpass filter (for example, as in the Luminex Guava easyCyte).