

**PerCP Anti-human CD144 Antibody  
\*55-7H1\***Catalog number: 114401S0, 114401S1, 114401S2  
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	PerCP

**Biological Properties**

---

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

**Spectral Properties**

---

Conjugate	PerCP
Excitation Wavelength	477 nm
Emission Wavelength	678 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 PerCP (ex/em = 477/678 nm). It is compatible with the 488 nm laser and 667/30 nm bandpass filter (for example, as in the Agilent

Technologies NovoCyte Advanteon).