

**APC Anti-human CD94 Antibody \*HP-3D9\***Catalog number: 109401C0, 109401C1, 109401C2  
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	CD94 (KLRD1, KP43)
Clone	HP-3D9
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**

---

The HP-3D9 monoclonal antibody binds with human CD94, a 43 kD glycoprotein typically located on the surface of T cells and natural killer cells. CD94 is associated with a variety of biologically interesting macromolecules/ligands, namely, NKG2-A, HLA class I and p39. CD94 is a fairly uncommon antibody target, with a little more than 2800 publications in the last decade. Even still, CD94 has a variety of applications in innate immunity and immunology research, commonly 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 702/87 nm bandpass filter (for example, as in the Luminex Amnis CellStream).