

PacOrange Anti-human CD94 Antibody
HP-3D9Catalog number: 109401L0, 109401L1
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	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	PacOrange

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PacOrange under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	PacOrange
Excitation Wavelength	400 nm
Emission Wavelength	551 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 PacOrange (ex/em = 400/551 nm). It is compatible with the 405 nm

laser and 525/45 nm bandpass filter (for example, as in the Agilent Technologies NovoCyte Advanteon).