

PE Anti-mouse CD26 Antibody *H194-112*Catalog number: 102601M0, 102601M1, 102601M2
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	Mouse
Class	Primary
Clonality	Monoclonal
Host	Rat
Isotype	Rat IgG2a kappa
Immunogen	CD26 (DPP IV ectoenzyme, ADA-binding protein, ADCP2)
Clone	H194-112
Conjugate	PE

Biological Properties

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

Spectral Properties

Conjugate	PE
Excitation Wavelength	566 nm
Emission Wavelength	574 nm

Applications

H194-112 is an anti-mouse monoclonal antibody that recognizes the CD26 antigen. CD26 (also known as DPP IV ectoenzyme, ADA-binding protein or ADCP2) is a 110 kD glycoprotein that is located on the surface of cells such as macrophages and NK cells. In some organisms, CD26 is an enhancer of cell population proliferation and plays a role in the downregulation of extracellular matrix disassembly, and is associated with a variety of biologically interesting macromolecules/ligands, namely, CD45, collagen and adenosine deaminase. CD26 is a fairly uncommon antibody target, with a little more than 3500 publications in the last decade. Even still, CD26 is frequently used in flow cytometry applications as a phenotypic marker for differentiation of cell types, particularly in the study of immunology. This antibody was purified through affinity chromatography and conjugated to PE (ex/em = 566/574 nm). It is compatible with the 561 nm laser and 586/20 nm bandpass filter (for

example, as in the Agilent Technologies NovoCyte).