

iFluor™ 350 Anti-mouse CD26 Antibody
H194-112Catalog number: 10260010, 10260011
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	Mouse
Class	Primary
Clonality	Monoclonal
Host	Rat
Isotype	Rat IgG2a kappa
Immunogen	CD26 (DPP IV ectoenzyme, ADA-binding protein, ADCP2)
Clone	H194-112
Conjugate	iFluor™ 350

Biological Properties

Appearance	Off-white liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 350 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 350
Excitation Wavelength	345 nm
Emission Wavelength	450 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 iFluor™ 350 (ex/em = 345/450 nm). It is compatible with the 355 nm laser and 450/50 nm bandpass filter (for example, as in the BD Special Order LSRFortessa™ Cell Analyzer).