

XFD594 Anti-mouse CD26 Antibody *H194-112*

Catalog Number: 10260170, 10260171

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	Lot specific (please consult certificate of analysis for given lot)
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	AF594

Biological Properties

Appearance	Purple liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with AF594 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging
Recommended Dilutions	For flow cytometry applications, the suggested concentration is at 5 uL/million cells in 100 uL staining buffer. For the best performance of each application, the optimal concentration of this reagent needs to be carefully determined. <i>*The suggested working dilution is provided as a guide only. It is recommended that the users titrates the product for use in their tests using proper positive and negative controls.</i>

Spectral Properties

Conjugate	AF594
Excitation Wavelength	590 nm
Emission Wavelength	618 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 XFD594 (ex/em = 590/618 nm). XFD594 is manufactured by AAT Bioquest, and it has a chemical structure similar to that of Alexa Fluor® 594 (Alexa Fluor® is the trademark of Thermo Fisher). It is compatible with the 592 nm laser and 610/30 nm bandpass filter (for example, as in the Luminex Amnis ImageStream).