

**XFD594 Anti-human CD209 Antibody \*UW60.1\***

Catalog Number: 12090170, 12090171

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	Human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Isotype	Mouse IgG1
Immunogen	CD209 (DC-SIGN)
Clone	UW60.1
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

---

UW60.1 is an anti-human monoclonal antibody that targets the CD209 antigen. CD209 (alternatively called DC-SIGN) is a transmembrane protein that is located on the surface of cells like macrophages, endothelial cells and dendritic cells. CD209 is associated with a variety of biologically interesting macromolecules/ligands, for example, mannose-bearing glycoproteins on several pathogens including HIV gp120. CD209 is a fairly uncommon antibody target, with a little more than 2700 publications in the last decade. Even still, CD209 is frequently used in flow cytometry applications as a phenotypic marker for differentiation of cell types, especially in the study of innate immunity and 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).