

iFluor™ A7 Anti-human CD209 Antibody
UW60.1Catalog number: 120900S0, 120900S1
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
Immunogen	CD209 (DC-SIGN)
Clone	UW60.1
Conjugate	iFluor™ A7

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ A7 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ A7
-----------	------------

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 iFluor™ A7.