

**iFluor™ 660 Anti-human CD209 Antibody**  
**\*UW60.1\***Catalog number: 120900G0, 120900G1  
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™ 660     |

**Biological Properties**

|             |                                                                                                            |
|-------------|------------------------------------------------------------------------------------------------------------|
| Appearance  | Blue liquid                                                                                                |
| Preparation | Antibody purified by affinity chromatography and then conjugated with iFluor™ 660 under optimal conditions |
| Application | Flow Cytometry (FACS), Fluorescence Imaging                                                                |

**Spectral Properties**

|                       |             |
|-----------------------|-------------|
| Conjugate             | iFluor™ 660 |
| Excitation Wavelength | 663 nm      |
| Emission Wavelength   | 678 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 iFluor™ 660 (ex/em = 663/678 nm). It is compatible with the 638 nm laser and 660/20 nm bandpass filter (for example, as in the Beckman Coulter Navios EX).