

**iFluor™ 710 Anti-human CD205 Antibody**  
**\*HD30\***Catalog number: 120500K0, 120500K1  
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 kappa |
| Immunogen          | CD205 (DEC205)   |
| Clone              | HD30             |
| Conjugate          | iFluor™ 710      |

**Biological Properties**

|             |                                                                                                            |
|-------------|------------------------------------------------------------------------------------------------------------|
| Appearance  | Blue liquid                                                                                                |
| Preparation | Antibody purified by affinity chromatography and then conjugated with iFluor™ 710 under optimal conditions |
| Application | Flow Cytometry (FACS), Fluorescence Imaging                                                                |

**Spectral Properties**

|                       |             |
|-----------------------|-------------|
| Conjugate             | iFluor™ 710 |
| Excitation Wavelength | 717 nm      |
| Emission Wavelength   | 739 nm      |

**Applications**

HD30 is an anti-human monoclonal antibody that targets the CD205 antigen. CD205 (alternatively called DEC205) is a 210 kD glycoprotein that is found on the surface of cells such as B cells. CD205 is associated with a variety of biologically interesting macromolecules/ligands, for example, MIR98. CD205 is a fairly uncommon antibody target, with a little more than 1200 publications in the last decade. Even still, CD205 has a variety

of applications in immunology research, often serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to iFluor™ 710 (ex/em = 717/739 nm).