

**iFluor™ 633 Anti-human CD53 Antibody**  
**\*HI29\***Catalog number: 105300E0, 105300E1  
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          | CD53 (Tetraspanin-25, MOX44) |
| Clone              | HI29                         |
| Conjugate          | iFluor™ 633                  |

**Biological Properties**

|             |                                                                                                            |
|-------------|------------------------------------------------------------------------------------------------------------|
| Appearance  | Black liquid                                                                                               |
| Preparation | Antibody purified by affinity chromatography and then conjugated with iFluor™ 633 under optimal conditions |
| Application | Flow Cytometry (FACS), Fluorescence Imaging                                                                |

**Spectral Properties**

|                       |             |
|-----------------------|-------------|
| Conjugate             | iFluor™ 633 |
| Excitation Wavelength | 640 nm      |
| Emission Wavelength   | 654 nm      |

**Applications**

HI29 is an anti-human monoclonal antibody that targets the CD53 antigen. CD53 (sometimes referred to as Tetraspanin-25, MOX44 or OX44) is a 35 - 42 kD member of the tetraspan family that is found on the surface of cells such as NK cells. In some organisms, CD53 enhances myoblast fusion, and is associated with a variety of biologically interesting macromolecules/ligands, in particular, VLA-4, Integrins and HLA-DR. CD53 is a

relatively rare antibody target, with fewer than 600 publications in the last decade. Even still, CD53 has a variety of applications in costimulatory molecules 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™ 633 (ex/em = 640/654 nm). It is compatible with the 640 nm laser and 660/20 nm bandpass filter (for example, as in the Agilent Technologies NovoCyte).