

**PE/iFluor™ 700 Anti-human CD93 Antibody**  
**\*VIMD2\***Catalog number: 109301X0, 109301X1, 109301X2  
Unit size: 25 tests, 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          | CD93 (C1QR1, MXRA4) |
| Clone              | VIMD2               |
| Conjugate          | PE/iFluor™ 700      |

**Biological Properties**

|             |                                                                                                               |
|-------------|---------------------------------------------------------------------------------------------------------------|
| Preparation | Antibody purified by affinity chromatography and then conjugated with PE/iFluor™ 700 under optimal conditions |
| Application | Flow Cytometry (FACS)                                                                                         |

**Spectral Properties**

|                       |                |
|-----------------------|----------------|
| Conjugate             | PE/iFluor™ 700 |
| Excitation Wavelength | 566 nm         |
| Emission Wavelength   | 708 nm         |

**Applications**

The VIMD2 monoclonal antibody reacts with human CD93, a 110 kD transmembrane glycoprotein often found on the surface of macrophages, monocytes, endothelial cells, platelets and granulocytes. CD93 is associated with a variety of biologically interesting macromolecules/ligands. CD93 is a fairly uncommon antibody target, with a little more than 1100 publications in the last decade. Even still, CD93 has a variety of applications in immunology research, frequently serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to PE/iFluor™ 700 (ex/em = 566/708 nm). It is compatible with the

561 nm laser and 710/50 nm bandpass filter (for example, as in the BD Special Order LSRFortessa™ Cell Analyzer).