

**FITC Anti-human CD87 Antibody \*VIM5\***Catalog number: 10870110, 10870111  
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          | CD87 (UPAR, PLAUR) |
| Clone              | VIM5               |
| Conjugate          | FITC               |

**Biological Properties**

|             |                                                                                                     |
|-------------|-----------------------------------------------------------------------------------------------------|
| Preparation | Antibody purified by affinity chromatography and then conjugated with FITC under optimal conditions |
| Application | Flow Cytometry (FACS), Fluorescence Imaging                                                         |

**Spectral Properties**

|                       |        |
|-----------------------|--------|
| Conjugate             | FITC   |
| Excitation Wavelength | 491 nm |
| Emission Wavelength   | 516 nm |

**Applications**

The VIM5 monoclonal antibody binds with human CD87, a 36 - 68 kD transmembrane protein often located on the surface of granulocytes and keratinocytes. CD87 is associated with a variety of biologically interesting macromolecules/ligands, in particular, Pro-UPA. CD87 is a relatively rare antibody target, with fewer than 400 publications in the last decade. Even still, CD87 has been widely used in immunology research, typically serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to FITC (ex/em = 491/516 nm). It is compatible with the 488 nm laser and 512/18 nm bandpass filter (for example, as in the Luminex Guava easyCyte).