

Purified Anti-human CD87 Antibody *VIM5*

Catalog number: 10870000

Unit size: 100 ug

Product Details

| | |
|--------------------|---|
| Storage Conditions | 2-8°C with minimized light exposure. Do not freeze. |
| Expiration Date | 12 months upon receiving |
| Concentration | 0.5 mg/mL |
| Formulation | Phosphate-buffered saline (PBS, pH 7.2), 0.09% sodium azide |

Antibody Properties

| | |
|--------------------|--------------------|
| Species Reactivity | Human |
| Class | Primary |
| Clonality | Monoclonal |
| Host | Mouse |
| Isotype | Mouse IgG1 |
| Immunogen | CD87 (UPAR, PLAUR) |
| Clone | VIM5 |

Biological Properties

| | |
|-------------|--|
| Appearance | Liquid |
| Preparation | Antibody purified by affinity chromatography |
| Application | Flow Cytometry (FACS), ELISA, HC, Western Blot |

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.