

**Purified Anti-human CD57 Antibody**  
**\*HI57a\***Catalog number: 10570000  
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 IgM           |
| Immunogen          | CD57 (HNK-1, Leu-7) |
| Clone              | HI57a               |

**Biological Properties**

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

**Applications**

The HI57a monoclonal antibody binds with human CD57, a 100 - 115 kD transmembrane protein often expressed on the surface of T cells, natural killer cells, granulocytes and platelets. CD57 is associated with a variety of biologically interesting macromolecules/ligands, for example, P-Selectin and Laminin. CD57 is a fairly uncommon antibody target, with a little more than 3000 publications in the last decade. Even still, CD57 has a variety of applications in costimulatory molecules and immunology research, frequently serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography.