

**Purified Anti-human CD300a Antibody**  
**\*MEM-260\***Catalog number: 13000000  
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          | CD300a (CLM-8, IRp60, CMRF-35H) |
| Clone              | MEM-260                         |

**Biological Properties**

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

**Applications**

MEM-260 is an anti-human monoclonal antibody that is specific for the CD300a antigen. CD300a (alternatively called CLM-8) is a transmembrane protein that is expressed on the surface of cells such as macrophages, T cells, dendritic cells and B cells. In some organisms, CD300a is a repressor of MyD88-dependent toll-like receptor signaling pathway, suppresses B cell receptor signaling pathway and is a promoter of phosphoprotein phosphatase activity. Also, it is a member of important cellular pathways, for example, the negative regulation of B cell receptor signaling pathway, negative regulation of MyD88-dependent toll-like receptor signaling pathway and regulation of T cell receptor signaling pathway. From a research standpoint, it is of biological interest due to its association with critical macromolecules/ligands like unknown. CD300a is a relatively rare antibody target, with fewer than 400 publications in the last decade. Even still, CD300a is essential for innate immunity and immunology research, commonly serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography.