

Purified Anti-human CD50 Antibody
MEM-04Catalog number: 10500000
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 IgG2b kappa |
| Immunogen | CD50 (ICAM-3) |
| Clone | MEM-04 |

Biological Properties

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

Applications

MEM-04 is an anti-human monoclonal antibody that targets the CD50 antigen. CD50 (alternatively called ICAM-3) is a 130 kD member of the Ig superfamily that is located on the surface of cells like macrophages. CD50 is associated with a variety of biologically interesting macromolecules/ligands, in particular, integrin α d/ β 2 and LFA-1. CD50 is a relatively rare antibody target, with fewer than 300 publications in the last decade. Even still, CD50 is typically used in flow cytometry applications as a phenotypic marker for differentiation of cell types, particularly in the study of costimulatory molecules and cell biology. This antibody was purified through affinity chromatography.