

**Purified Anti-dog/ chicken/ rabbit/ guinea pig/ horse/ cow/ mouse/ rat/ pig/ non-human primates/ human CD79a Antibody \*HM47\***

Catalog number: 10790000  
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 | Dog, chicken, rabbit, guinea pig, horse, cow, mouse, rat, pig, non-human primates, human |
| Class              | Primary                                                                                  |
| Clonality          | Monoclonal                                                                               |
| Host               | Mouse                                                                                    |
| Isotype            | Mouse IgG1 kappa                                                                         |
| Immunogen          | CD79a (Mb-1, IGA)                                                                        |
| Clone              | HM47                                                                                     |

### Biological Properties

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

### Applications

HM47 is an anti-dog/ chicken/ rabbit/ guinea pig/ horse/ cow/ mouse/ rat/ pig/ non-human primates/ human monoclonal antibody that forms an immune complex with the CD79a antigen. CD79a (sometimes referred to as Mb-1 or IGA) is a 47 kD single-pass type I membrane protein that is expressed on the surface of cells such as B cells. CD79a is a component of essential cellular pathways, in particular, the B cell receptor signaling pathway. From a research standpoint, it is of biological interest due to its association with key macromolecules/ligands like CD22, CD79b, CD5 and CD19. CD79a is a fairly uncommon antibody target, with a little more than 4100 publications in the last decade. Even still, CD79a has been widely used in immunology research, commonly serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography.