

**mFluor™ UV460 Anti-dog/ chicken/ rabbit/  
guinea pig/ horse/ cow/ mouse/ rat/ pig/  
non-human primates/ human CD79a  
Antibody \*HM47\***

Catalog number: 107900Y0, 107900Y1  
Unit size: 100 tests, 500 tests

## Product Details

|                    |                                                                             |
|--------------------|-----------------------------------------------------------------------------|
| Storage Conditions | 2-8°C with minimized light exposure. Do not freeze.                         |
| Expiration Date    | 12 months upon receiving                                                    |
| Concentration      | 0.1 mg/mL                                                                   |
| Formulation        | Phosphate-buffered saline (PBS, pH 7.2), 0.09% sodium azide, 0.2% (w/v) BSA |

## 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                                                                                     |
| Conjugate          | mFluor™ UV460                                                                            |

## Biological Properties

|             |                                                                                                              |
|-------------|--------------------------------------------------------------------------------------------------------------|
| Preparation | Antibody purified by affinity chromatography and then conjugated with mFluor™ UV460 under optimal conditions |
| Application | Flow Cytometry (FACS), Fluorescence Imaging                                                                  |

## Spectral Properties

|                       |               |
|-----------------------|---------------|
| Conjugate             | mFluor™ UV460 |
| Excitation Wavelength | 358 nm        |
| Emission Wavelength   | 456 nm        |

## 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 and conjugated to mFluor™ UV460 (ex/em = 358/456 nm). It is compatible with the 355 nm laser and 450/50 nm bandpass filter (for example, as in the BD LSRFortessa™ Cell Analyzer).