

**iFluor™ 820 Anti-dog/ chicken/ rabbit/
guinea pig/ horse/ cow/ mouse/ rat/ pig/
non-human primates/ human CD79a
Antibody *HM47***

Catalog number: 107900P0, 107900P1
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	iFluor™ 820

Biological Properties

Appearance	Green liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 820 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 820
Excitation Wavelength	822 nm
Emission Wavelength	850 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 iFluor™ 820 (ex/em = 822/850 nm).