

PE Anti-human CD93 Antibody *VIMD2*Catalog number: 109301M0, 109301M1, 109301M2
Unit size: 25 tests, 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	Human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Isotype	Mouse IgG1
Immunogen	CD93 (C1QR1, MXRA4)
Clone	VIMD2
Conjugate	PE

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PE under optimal conditions
Application	Flow Cytometry (FACS)

Spectral Properties

Conjugate	PE
Excitation Wavelength	566 nm
Emission Wavelength	574 nm

Applications

The VIMD2 monoclonal antibody reacts with human CD93, a 110 kD transmembrane glycoprotein often found on the surface of macrophages, monocytes, endothelial cells, platelets and granulocytes. CD93 is associated with a variety of biologically interesting macromolecules/ligands. CD93 is a fairly uncommon antibody target, with a little more than 1100 publications in the last decade. Even still, CD93 has a variety of applications in immunology research, frequently serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to PE (ex/em = 566/574 nm). It is compatible with the 561 nm laser and 586/20 nm bandpass filter (for example, as in the Agilent Technologies NovoCytte Quanteon).