

mFluor™ UV375 Anti-human CD93 Antibody
VIMD2Catalog number: 109300X0, 109300X1
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	Human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Isotype	Mouse IgG1
Immunogen	CD93 (C1QR1, MXRA4)
Clone	VIMD2
Conjugate	mFluor™ UV375

Biological Properties

Appearance	Yellow liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with mFluor™ UV375 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	mFluor™ UV375
Excitation Wavelength	351 nm
Emission Wavelength	387 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 mFluor™ UV375 (ex/em = 351/387 nm). It is compatible with the 355 nm laser and 379/28 nm bandpass filter (for example, as in the BD FACSymphony™ A5).