

iFluor™ 546 Anti-human CD87 Antibody
VIM5Catalog number: 10870080, 10870081
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	CD87 (UPAR, PLAUR)
Clone	VIM5
Conjugate	iFluor™ 546

Biological Properties

Appearance	Red liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 546 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 546
Excitation Wavelength	541 nm
Emission Wavelength	557 nm

Applications

The VIM5 monoclonal antibody binds with human CD87, a 36 - 68 kD transmembrane protein often located on the surface of granulocytes and keratinocytes. CD87 is associated with a variety of biologically interesting macromolecules/ligands, in particular, Pro-UPA. CD87 is a relatively rare antibody target, with fewer than 400 publications in the last decade. Even still, CD87 has been widely used in immunology research,

typically serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to iFluor™ 546 (ex/em = 541/557 nm). It is compatible with the 532 nm laser and 575/25 nm bandpass filter (for example, as in the BD Special Order LSRFortessa™ Cell Analyzer).