

iFluor™ 555 Anti-human CD109 Antibody
W7C5Catalog number: 11090090, 11090091
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	CD109 (CPAMD7, Gov platelet alloantigen)
Clone	W7C5
Conjugate	iFluor™ 555

Biological Properties

Appearance	Red liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 555 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 555
Excitation Wavelength	557 nm
Emission Wavelength	570 nm

Applications

The W7C5 monoclonal antibody binds with human CD109, a 162 kD transmembrane protein typically expressed on the surface of T cells and T cell lines. CD109 acts in essential cellular pathways, for instance, the negative regulation of transforming growth factor beta receptor signaling pathway. Additionally, in some organisms, it is a suppressor of transforming growth factor beta receptor signaling pathway, is a repressor of

wound healing and is a repressor of keratinocyte proliferation. From a research standpoint, it is of biological interest due to its association with critical macromolecules/ligands. CD109 is a relatively rare antibody target, with fewer than 600 publications in the last decade. Even still, CD109 is vital to 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 iFluor™ 555 (ex/em = 557/570 nm). It is compatible with the 561 nm laser and 577/35 nm bandpass filter (for example, as in the Luminex Amnis ImageStream).