

**PE/iFluor™ 594 Anti-human CD109 Antibody
*W7C5***Catalog number: 110901Y0, 110901Y1, 110901Y2
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 | CD109 (CPAMD7, Gov platelet alloantigen) |
| Clone | W7C5 |
| Conjugate | PE/iFluor™ 594 |

Biological Properties

| | |
|-------------|---|
| Preparation | Antibody purified by affinity chromatography and then conjugated with PE/iFluor™ 594 under optimal conditions |
| Application | Flow Cytometry (FACS) |

Spectral Properties

| | |
|-----------------------|----------------|
| Conjugate | PE/iFluor™ 594 |
| Excitation Wavelength | 566 nm |
| Emission Wavelength | 606 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 PE/iFluor™ 594 (ex/em = 566/606 nm). It is compatible with the 561 nm laser and 610/30 nm bandpass filter (for example, as in the Luminex Amnis FlowSight).