

**iFluor™ 840 Anti-human CD9 Antibody
*HI9a***Catalog number: 100900Q0, 100900Q1
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 | CD9 (MRP-1, P24) |
| Clone | HI9a |
| Conjugate | iFluor™ 840 |

Biological Properties

| | |
|-------------|--|
| Appearance | Dark brown liquid |
| Preparation | Antibody purified by affinity chromatography and then conjugated with iFluor™ 840 under optimal conditions |
| Application | Flow Cytometry (FACS), Fluorescence Imaging |

Spectral Properties

| | |
|-----------------------|-------------|
| Conjugate | iFluor™ 840 |
| Excitation Wavelength | 836 nm |
| Emission Wavelength | 879 nm |

Applications

HI9a is an anti-human monoclonal antibody that recognizes the CD9 antigen. CD9 (alternatively called MIC3, TSPAN29, Tetraspanin or 5H9 antigen) is a 24 kD member of the Tetraspan family that is expressed on the surface of cells like macrophages, granulocytes, epithelial cells, platelets and endothelial cells. In many organisms, CD9 is involved in the negative regulation of cellular component movement, represses cell

population proliferation and negatively regulates platelet aggregation. From a research standpoint, it is of biological interest due to its association with critical macromolecules/ligands such as Integrin β 1. CD9 is a fairly uncommon antibody target, with a little more than 6100 publications in the last decade. Even still, CD9 is frequently used in flow cytometry applications as a phenotypic marker for differentiation of cell types, particularly in the study of immunology and stem cells. This antibody was purified through affinity chromatography and conjugated to iFluor™ 840 (ex/em = 836/879 nm).