

**iFluor™ A7 Anti-human CD41 Antibody**  
**\*HIP8\***Catalog number: 104100S0, 104100S1  
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          | CD41 (GPIIb, ITGA2B) |
| Clone              | HIP8                 |
| Conjugate          | iFluor™ A7           |

**Biological Properties**

|             |                                                                                                           |
|-------------|-----------------------------------------------------------------------------------------------------------|
| Preparation | Antibody purified by affinity chromatography and then conjugated with iFluor™ A7 under optimal conditions |
| Application | Flow Cytometry (FACS), Fluorescence Imaging                                                               |

**Spectral Properties**

|           |            |
|-----------|------------|
| Conjugate | iFluor™ A7 |
|-----------|------------|

**Applications**

HIP8 is an anti-human monoclonal antibody that recognizes the CD41 antigen. CD41 (also known as GP2B) is a 22 kD member of the Integrin family that is found on the surface of cells such as stem cells. CD41 is a component of critical cellular pathways, for instance, the integrin-mediated signaling pathway. Moreover, in some organisms, it enhances leukocyte migration. From a research standpoint, it is of biological interest due to its association with key macromolecules/ligands such as von Willebrand factor (vWF) and Fibrinogen. CD41 is a fairly uncommon antibody target, with a little more than 4000 publications in the last decade. Even still, CD41 is frequently used in flow cytometry applications as a phenotypic marker for differentiation of cell types, specifically in the study of cell biology and cell adhesion. This antibody was purified through affinity chromatography and conjugated to iFluor™ A7.