

**mFluor™ UV375 Anti-human CD150  
Antibody \*SLAM.4\***Catalog number: 115000X0, 115000X1  
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          | CD150 (SLAM, IPO-3) |
| Clone              | SLAM.4              |
| Conjugate          | mFluor™ UV375       |

**Biological Properties**

|             |                                                                                                              |
|-------------|--------------------------------------------------------------------------------------------------------------|
| Appearance  | Yellow liquid                                                                                                |
| Preparation | Antibody purified by affinity chromatography and then conjugated with mFluor™ UV375 under optimal conditions |
| Application | Flow Cytometry (FACS), Fluorescence Imaging                                                                  |

**Spectral Properties**

|                       |               |
|-----------------------|---------------|
| Conjugate             | mFluor™ UV375 |
| Excitation Wavelength | 351 nm        |
| Emission Wavelength   | 387 nm        |

**Applications**

The SLAM.4 monoclonal antibody binds with human CD150, a 70 - 95 kD glycoprotein commonly located on the surface of B cells, dendritic cells, endothelial cells, T cells and Tregs. CD150 plays a role in important cellular pathways, in particular, the negative regulation of CD40 signaling pathway. In addition, in certain organisms, it is a negative regulator of tumor necrosis factor production, suppresses interleukin-6 production

and plays a role in the upregulation of JNK cascade. From a research standpoint, it is of biological interest due to its association with critical macromolecules/ligands such as tyrosine phosphatase CD45. CD150 is a fairly uncommon antibody target, with a little more than 2700 publications in the last decade. Even still, CD150 is often used in flow cytometry applications as a phenotypic marker for differentiation of cell types, particularly in the study of immunology. This antibody was purified through affinity chromatography and conjugated to mFluor™ UV375 (ex/em = 351/387 nm). It is compatible with the 355 nm laser and 379/28 nm bandpass filter (for example, as in the BD LSRFortessa™ X-20).