

**iFluor™ 560 Anti-human CD107 Antibody**  
**\*H4B4\***Catalog number: 110710A0, 110710A1  
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 kappa      |
| Immunogen          | CD107b (LAMP2, LAMPb) |
| Clone              | H4B4                  |
| Conjugate          | iFluor™ 560           |

**Biological Properties**

|             |                                                                                                            |
|-------------|------------------------------------------------------------------------------------------------------------|
| Preparation | Antibody purified by affinity chromatography and then conjugated with iFluor™ 560 under optimal conditions |
| Application | Flow Cytometry (FACS), Fluorescence Imaging                                                                |

**Spectral Properties**

|                       |             |
|-----------------------|-------------|
| Conjugate             | iFluor™ 560 |
| Excitation Wavelength | 560 nm      |
| Emission Wavelength   | 571 nm      |

**Applications**

H4B4 is an anti-human monoclonal antibody that targets the CD107b antigen. CD107b (sometimes called LAMP2 or LAMPb) is a 45 kD transmembrane protein that is found on the surface of cells like granulocytes and endothelial cells. In many organisms, CD107 is a suppressor of protein-containing complex assembly. Additionally, it has been associated with essential biological processes like chaperone-mediated autophagy, particularly protein targeting to lysosome involved in chaperone-mediated autophagy. From a research standpoint, it is of biological interest due to its association with vital macromolecules/ligands. CD107 is a relatively rare antibody target, with fewer than 800 publications in

the last decade. Even still, CD107b is typically used in flow cytometry applications as a phenotypic marker for differentiation of cell types, specifically in the study of protein trafficking and clearance and neuroscience. This antibody was purified through affinity chromatography and conjugated to iFluor™ 560 (ex/em = 560/571 nm). It is compatible with the 561 nm laser and 586/15 nm bandpass filter (for example, as in the BD FACSCelesta™).