

**PE/iFluor™ 700 Anti-human CD107 Antibody
*H4B4***Catalog number: 110711U0, 110711U1, 110711U2
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 kappa
Immunogen	CD107b (LAMP2, LAMPb)
Clone	H4B4
Conjugate	PE/iFluor™ 700

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PE/iFluor™ 700 under optimal conditions
Application	Flow Cytometry (FACS)

Spectral Properties

Conjugate	PE/iFluor™ 700
Excitation Wavelength	566 nm
Emission Wavelength	708 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 PE/iFluor™ 700 (ex/em = 566/708 nm). It is compatible with the 561 nm laser and 710/50 nm bandpass filter (for example, as in the BD LSRFortessa™ X-20).