

**iFluor™ 660 Anti-human/ non-human
primates/ mouse CD107a Antibody *H4A3***Catalog number: 110700G0, 110700G1
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, non-human primates, mouse
Class	Primary
Clonality	Monoclonal
Host	Mouse
Isotype	Mouse IgG1 kappa
Immunogen	CD107a (LAMP-1, LGP-120)
Clone	H4A3
Conjugate	iFluor™ 660

Biological Properties

Appearance	Blue liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 660 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 660
Excitation Wavelength	663 nm
Emission Wavelength	678 nm

Applications

The H4A3 monoclonal antibody binds with human/ non-human primates/ mouse CD107a, a 45 kD transmembrane protein frequently expressed on the surface of macrophages, endothelial cells, epithelial cells and activated platelets. CD107a is associated with a variety of biologically interesting macromolecules/ligands, for example, . CD107a is a fairly uncommon antibody target, with a little more than 4600 publications in

the last decade. Even still, CD107a is often used in flow cytometry applications as a phenotypic marker for differentiation of cell types, particularly in the study of immunology and neuroscience. This antibody was purified through affinity chromatography and conjugated to iFluor™ 660 (ex/em = 663/678 nm). It is compatible with the 640 nm laser and 667/30 nm bandpass filter (for example, as in the Agilent Technologies NovoCyt^e Advanteon).