

iFluor™ A7 Anti-human/ non-human primates/ mouse CD107a Antibody *H4A3*Catalog number: 110700S0, 110700S1
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™ 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

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™ A7.