

PE/iFluor™ 750 Anti-human CD98 Antibody
MEM-108Catalog number: 109801T0, 109801T1, 109801T2
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
Immunogen	CD98 (SLC3A2, MDU1, 4F2hc, RL-388, FRP-1, 4F2)
Clone	MEM-108
Conjugate	PE/iFluor™ 750

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PE/iFluor™ 750 under optimal conditions
Application	Flow Cytometry (FACS)

Spectral Properties

Conjugate	PE/iFluor™ 750
Excitation Wavelength	566 nm
Emission Wavelength	778 nm

Applications

MEM-108 is an anti-human monoclonal antibody that targets the CD98 antigen. CD98 (sometimes called 4F2hc, SLC3A2, MDU1 or RL-388) is a 80 kD transmembrane protein that is located on the surface of cells like platelets, endothelial cells, NK cells, B cells and epithelial cells. CD98 is associated with a variety of biologically interesting macromolecules/ligands, for instance, actin. CD98 is a fairly uncommon antibody target, with a little more than 1300 publications in the last decade. Even still, CD98 is often used in flow cytometry applications as a phenotypic marker for differentiation of cell types, specifically in the study of immunology. This antibody was purified through affinity chromatography and conjugated

to PE/iFluor™ 750 (ex/em = 566/778 nm). It is compatible with the 561 nm laser and 780/60 nm bandpass filter (for example, as in the BD FACS Aria™ III).