

TRITC Anti-human CD98 Antibody
MEM-108Catalog number: 109801J0, 109801J1
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
Immunogen	CD98 (SLC3A2, MDU1, 4F2hc, RL-388, FRP-1, 4F2)
Clone	MEM-108
Conjugate	TRITC

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with TRITC under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	TRITC
Excitation Wavelength	544 nm
Emission Wavelength	570 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 TRITC (ex/em = 544/570 nm). It is compatible with the 561 nm laser and 577/15 nm bandpass filter (for example, as in the Bio-Rad ZE5 Cell Analyzer).