

iFluor™ 430 Anti-human CD98 Antibody
MEM-108Catalog number: 10980030, 10980031
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	iFluor™ 430

Biological Properties

Appearance	Yellow liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 430 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 430
Excitation Wavelength	433 nm
Emission Wavelength	498 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 iFluor™ 430 (ex/em = 433/498 nm). It is compatible with the 445 nm laser and 510/80 nm bandpass filter (for example, as in the BD FACSAria™ Fusion).