

iFluor™ 800 Anti-human CD97 Antibody
MEM-180Catalog number: 109700N0, 109700N1
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	CD97 (BL-KDDF12)
Clone	MEM-180
Conjugate	iFluor™ 800

Biological Properties

Appearance	Green liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 800 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 800
Excitation Wavelength	801 nm
Emission Wavelength	820 nm

Applications

The MEM-180 monoclonal antibody binds with human CD97, a 74 kD multi-pass membrane protein typically expressed on the surface of macrophages, dendritic cells, granulocytes and monocytes. CD97 is involved with essential cellular pathways, namely, the G protein-coupled receptor signaling pathway and cell surface receptor signaling pathway. From a research standpoint, it is of biological interest due to its

association with important macromolecules/ligands such as CD55 (DAF). CD97 is a relatively rare antibody target, with fewer than 600 publications in the last decade. Even still, CD97 is typically used in flow cytometry applications as a phenotypic marker for differentiation of cell types, particularly in the study of immunology. This antibody was purified through affinity chromatography and conjugated to iFluor™ 800 (ex/em = 801/820 nm).