

**mFluor™ Green 620 Anti-human CD97  
Antibody \*MEM-180\***Catalog number: 109700U0, 109700U1  
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	mFluor™ Green 620

**Biological Properties**

Appearance	Purple liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with mFluor™ Green 620 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

**Spectral Properties**

Conjugate	mFluor™ Green 620
Excitation Wavelength	525 nm
Emission Wavelength	623 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 mFluor™ Green 620 (ex/em = 525/623 nm). It is compatible with the 532 nm laser and 620/15 nm bandpass filter (for example, as in the Thermo Fisher Attune NxT).