

iFluor™ 790 Anti-human CD300a Antibody
MEM-260Catalog number: 130000M0, 130000M1
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	CD300a (CLM-8, IRp60, CMRF-35H)
Clone	MEM-260
Conjugate	iFluor™ 790

Biological Properties

Appearance	Green liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with iFluor™ 790 under optimal conditions
Application	Flow Cytometry (FACS), Fluorescence Imaging

Spectral Properties

Conjugate	iFluor™ 790
Excitation Wavelength	787 nm
Emission Wavelength	812 nm

Applications

MEM-260 is an anti-human monoclonal antibody that is specific for the CD300a antigen. CD300a (alternatively called CLM-8) is a transmembrane protein that is expressed on the surface of cells such as macrophages, T cells, dendritic cells and B cells. In some organisms, CD300a is a repressor of MyD88-dependent toll-like receptor signaling pathway, suppresses B cell receptor signaling pathway and is a promoter

of phosphoprotein phosphatase activity. Also, it is a member of important cellular pathways, for example, the negative regulation of B cell receptor signaling pathway, negative regulation of MyD88-dependent toll-like receptor signaling pathway and regulation of T cell receptor signaling pathway. From a research standpoint, it is of biological interest due to its association with critical macromolecules/ligands like unknown. CD300a is a relatively rare antibody target, with fewer than 400 publications in the last decade. Even still, CD300a is essential for innate immunity and immunology research, commonly serving as a phenotypic marker for differentiating cell types in flow cytometric applications. This antibody was purified through affinity chromatography and conjugated to iFluor™ 790 (ex/em = 787/812 nm).