

**PE Mouse Anti-human HLA-DR1 (empty)
Antibody *MEM-267, monoclonal***

Catalog number: V1031045

Unit size: 0.1 mg

Product Details

Storage Conditions	2-8°C with minimized light exposure. Do not freeze.
Expiration Date	12 months upon receiving
Concentration	Lot specific (please consult certificate of analysis for given lot)
Formulation	Phosphate-buffered saline (PBS, pH 7.2), 15 mM sodium azide, 0.2% (w/v) BSA

Antibody Properties

Species Reactivity	Human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	HLA-DR1 (empty)
Clone	MEM-267
Conjugate	PE

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PE under optimal conditions
Application	FC (QC TESTED)

Applications

HLA class II histocompatibility antigen, DR α chain is a 33 kDa transmembrane protein that can be found in the trans-Golgi network membrane, late endosome membrane and MHC class II protein complex of cells. In Homo sapiens, HLA class II histocompatibility antigen, DR α chain binds with peptide antigen, polysaccharide and MHC class II protein complex. It aids in organismal processes, namely, peptide antigen assembly with MHC class II protein complex, cognition and immune response. It has been closely linked to important functions like MHC class II receptor activity. Additionally, it is the subject of extensive examination due to the fact that it plays a role in the T cell receptor signaling pathway and interferon- γ -mediated signaling pathway. Sequencing of HLA class II histocompatibility antigen, DR α chain has supported it contains 3 conserved structural units: cytoplasmic, Ig-like C1-type and extracellular domain.