

Purified Azide Free Rat Anti-mouse MHC Class II Antibody *M5/114, monoclonal*

Catalog number: V1031760

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	Azide free Phosphate-buffered saline (PBS, pH 7.2) 0.2 µm filter sterilized, 0.2% (w/v) BSA

Antibody Properties

Species Reactivity	Mouse
Class	Primary
Clonality	Monoclonal
Host	Rat
Immunogen	MHC Class II
Clone	M5/114
Conjugate	Purified Azide Free

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with Purified Azide Free under optimal conditions
Application	FC (QC TESTED), IP, WB, IHC(P), IHC(F), FUNC

Applications

HLA class II histocompatibility antigen, DP α 1 chain is a 34 kDa transmembrane protein that can be found in the integral component of plasma membrane, transport vesicle membrane and ER to Golgi transport vesicle membrane of cells. It is sometimes called MHC class II DP3- α , MHC class II DPA1 and DP(W3). In Homo sapiens, HLA class II histocompatibility antigen, DP α 1 chain has been thought to be involved with vital functions like MHC class II receptor activity. It is an integral part of organismal processes, namely, adaptive immune response, immune response and cellular response to interferon- γ . HLA class II histocompatibility antigen, DP α 1 chain is the subject of extensive application because of the fact that it acts in the T cell receptor signaling pathway and interferon- γ -mediated signaling pathway. It enhances T cell proliferation, interferon- γ production and T cell activation, and moreover, binds to peptide antigen. Sequencing of HLA class II histocompatibility antigen, DP α 1 chain has supported it contains 3 conserved structural units: extracellular, cytoplasmic and Ig-like C1-type domain.