

PE Mouse Anti-human HLA-A2 Antibody
BB7.2, monoclonalCatalog number: V103885
Unit size: 100 tests**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-A2
Clone	BB7.2
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 I histocompatibility antigen, A α chain is a 47 kDa transmembrane protein that can be expressed in the integral component of luminal side of endoplasmic reticulum membrane, cell surface and recycling endosome membrane of cells. In Homo sapiens, HLA class I histocompatibility antigen, A α chain has been found to be involved in organismal processes, namely, T cell mediated cytotoxicity directed against tumor cell target, defense response to Gram-positive bacterium and CD8-positive, α -beta T cell activation. Sequencing of HLA class I histocompatibility antigen, A α chain has exemplified it contains 3 conserved structural units: extracellular, Ig-like C1-type and cytoplasmic domain. It is the subject of intensive research due to the fact that it plays a role in the antigen processing and presentation of endogenous peptide antigen via MHC class I via ER pathway, TAP-independent, TAP-dependent and type I interferon signaling pathway. Moreover, HLA class I histocompatibility antigen, A α chain recognizes signaling receptor, TAP complex and T cell receptor. It is a positive regulator of interferon- γ production, T cell mediated cytotoxicity and CD8-, α -beta T cell proliferation. It takes part in processes such as immune response.