

PE Mouse Anti-human/non-human primates HLA-B7 Antibody *BB7.1, monoclonal*Catalog number: V103905
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, non-human primates
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	HLA-B7
Clone	BB7.1
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, B α chain is a 47 kDa transmembrane protein that can be expressed in the early endosome membrane, recycling endosome membrane and extracellular exosome of cells. In humans, HLA class I histocompatibility antigen, B α chain takes part in processes such as interleukin-12 production, immune response and dendritic cell differentiation. It acts to positively regulate T cell mediated cytotoxicity. It reacts with chaperone, peptide antigen and TAP. Sequencing of HLA class I histocompatibility antigen, B α chain has exemplified it contains 3 conserved structural units: cytoplasmic, Ig-like C1-type and extracellular domain. HLA class I histocompatibility antigen, B α chain is the subject of extensive examination stemming from the fact that it is a component of the type I interferon signaling pathway, interferon- γ -mediated signaling pathway and antigen processing and presentation of endogenous peptide antigen via MHC class I, TAP-independent, and moreover, takes part in processes such as organismal processes, for example, viral process, adaptive immune response and defense response. HLA class I histocompatibility antigen, B α chain is clinically significant because abnormalities in its function have been thought to be involved with diseases like Stevens-Johnson syndrome (SJS) and Spondyloarthritis 1 (SPDA1).