

**Purified Mouse Anti-human/pig HLA-DQ1/DQ3 Antibody \*HL-37, monoclonal\***

Catalog number: V103910

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, pig
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	HLA-DQ1/DQ3
Clone	HL-37

**Biological Properties**

---

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

**Applications**

---

HLA class II histocompatibility antigen, DQ  $\alpha$  1 chain (also known as DC-1  $\alpha$  chain, HLA-DCA, DC- $\alpha$  or MHC class II DQA1) is a transmembrane protein with a molecular weight of 32 kDa, located in the transport vesicle membrane, endosome membrane and membrane of cells. In humans, DC-1  $\alpha$  chain reacts with peptide antigen. DC-1  $\alpha$  chain has been closely linked to important functions like MHC class II receptor activity. It is the subject of intensive examination stemming from the fact that it is a component of the T cell receptor signaling pathway and interferon- $\gamma$ -mediated signaling pathway. Sequencing of DC-1  $\alpha$  chain has supported it contains 3 conserved structural units: Ig-like C1-type, cytoplasmic and extracellular domain. HLA class II histocompatibility antigen, DQ  $\alpha$  1 chain has been found to be involved in organismal processes, for example, immune response, adaptive immune response and antigen processing and presentation of exogenous peptide antigen via MHC class II.