

**Purified Low Endotoxin Mouse Anti-non-human primates/rat TCR  $\alpha/\beta$  Antibody  
\*R73, monoclonal\***

Catalog number: V1032225

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	Low endotoxin azide free Phosphate-buffered saline (PBS, pH 7.2), 0.2 $\mu$ m filter sterilized. Endotoxin level is less than 0.01 EU/ $\mu$ g of the protein, as determined by the LAL test., 0.2% (w/v) BSA

### Antibody Properties

---

Species Reactivity	Non-human primates, rat
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	TCR $\alpha/\beta$
Clone	R73
Conjugate	Purified Low Endotoxin

### Biological Properties

---

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

### Applications

---

T cell receptor  $\alpha$  chain constant is a 18 kDa transmembrane protein that can be expressed in the integral component of membrane, plasma membrane and T cell receptor complex of cells. In Homo sapiens, T cell receptor  $\alpha$  chain constant is the subject of intensive study due to the fact that it is a component of the T cell receptor signaling pathway. Sequencing of T cell receptor  $\alpha$  chain constant has shown it contains 2 conserved structural units: Ig-like C1-type and cytoplasmic domain, and furthermore, aids in organismal processes, in particular, adaptive immune response. T cell receptor  $\alpha$  chain constant is thought to be essential to immune response. Mutations and abnormalities in T cell receptor  $\alpha$  chain constant have been thought to be involved with a number of diseases, for instance, immunodeficiency 7 (IMD7). Immunodeficiency 7, an autosomal recessive inherited disorder characterized by recurrent infections, lymphadenopathy and autoimmunity, has especially been of interest to scientists.