

## Purified Low Endotoxin Mouse Anti-human TCR V $\beta$ 5.3-related Antibody \*MEM-262, monoclonal\*

Catalog number: V1032315

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	Human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	TCR V $\beta$ 5.3-related
Clone	MEM-262
Conjugate	Purified Low Endotoxin

### Biological Properties

---

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

### Applications

---

M1-specific T cell receptor  $\beta$  chain is a transmembrane protein with a molecular weight of 40 kDa, located in the integral component of membrane and plasma membrane of cells. In Homo sapiens, M1-specific T cell receptor  $\beta$  chain is an integral part of organismal processes, for example, adaptive immune response. Sequencing of M1-specific T cell receptor  $\beta$  chain has demonstrated that it contains 3 conserved structural units: Ig-like C1-type, Ig-like V-type and cytoplasmic domain.