

**Purified Mouse Anti-human TCR  
Vβ5.3-related Antibody \*MEM-262,  
monoclonal\***

Catalog number: V1032310

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
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	TCR Vβ5.3-related
Clone	MEM-262

**Biological Properties**

---

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

**Applications**

---

M1-specific T cell receptor β 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 β chain is an integral part of organismal processes, for example, adaptive immune response. Sequencing of M1-specific T cell receptor β chain has demonstrated that it contains 3 conserved structural units: Ig-like C1-type, Ig-like V-type and cytoplasmic domain.