

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

Catalog number: V1032305

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	Azide free Phosphate-buffered saline (PBS, pH 7.2) 0.2 µm filter sterilized, 0.2% (w/v) BSA

Antibody Properties

Species Reactivity	Human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	TCR Vβ5.3-related
Clone	MEM-262
Conjugate	Purified Azide Free

Biological Properties

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

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.