

Biotin Mouse Anti-human TCR α/β Antibody
IP26, monoclonalCatalog number: V1032235
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 α/β
Clone	IP26
Conjugate	Biotin

Biological Properties

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

Applications

T cell receptor α 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 α 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 α 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 α chain constant is thought to be essential to immune response. Mutations and abnormalities in T cell receptor α chain constant have been thought to be involved with a number of diseases, for instance, immunodeficiency 7 (IMD7). Immunodeficiency 7, an autosomal recessive inheritance disorder characterized by recurrent infections, lymphadenopathy and autoimmunity, has especially been of interest to scientists.