

**Purified Mouse Anti-human IL-9 Antibody**  
**\*MH9A4, monoclonal\***Catalog number: V1031560  
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	IL-9
Clone	MH9A4

**Biological Properties**

Preparation	Antibody purified by affinity chromatography and then conjugated with under optimal conditions
Application	ELISA

**Applications**

Interleukin-9 (IL-9) is a 16 kDa protein that can be found in the extracellular region and extracellular space of cells. It is alternatively called T-cell growth factor P40 and Cytokine P40. In humans, IL-9 has been thought to be involved with critical functions like growth factor and cytokine activity. IL-9 reacts with interleukin-9 receptor. It is the subject of comprehensive study due to the fact that it acts in the interleukin-9-mediated signaling pathway. IL-9 is thought to be essential to organismal processes, for instance, inflammatory response and immune response. It plays a role in the upregulation of interleukin-5 biosynthetic process, cell population proliferation and cell growth.