

**PE Mouse Anti-human/mouse FoxP3  
Antibody \*3G3, monoclonal\***Catalog number: V103665  
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, mouse
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	FoxP3
Clone	3G3
Conjugate	PE

**Biological Properties**

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

**Applications**

Forkhead box protein P3 is a 47 kDa protein that can be found in the nuclear chromatin, cytoplasm and nucleoplasm of cells. In Homo sapiens, forkhead box protein P3 is the subject of intensive application due to the fact that it is a component of the T cell receptor signaling pathway and regulation of Wnt signaling pathway. Forkhead box protein P3 is a promoter of peripheral T cell tolerance induction, histone acetylation and DNA-templated transcription, on the other hand, also negatively regulates interleukin-17 production, interleukin-4 production and activated T cell proliferation. It is an integral part of DNA-templated transcription, isotype switching to IgG isotypes and T cell anergy, and also, binds to Forkhead and sequence-specific DNA. Forkhead box protein P3 has been associated with key functions like protein homodimerization, DNA-binding transcription factor and transcription corepressor activity, and furthermore, plays an important role in organismal processes, for instance, T cell mediated immunity, T cell activation and response to virus. Forkhead box protein P3 is clinically significant because abnormalities in its function have been thought to be involved with diseases such as Immunodeficiency polyendocrinopathy, enteropathy, X-linked syndrome (IPEX). Immunodeficiency polyendocrinopathy, enteropathy, X-linked syndrome, a x-linked recessive inherited disorder characterized by Diabetes mellitus, Immune dysregulation and Eosinophilia- has especially been of interest to researchers.