

**Purified Mouse Anti-human hGH Antibody**  
**\*GH-45, monoclonal\***Catalog number: V103815  
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	hGH
Clone	GH-45

**Biological Properties**

Preparation	Antibody purified by affinity chromatography and then conjugated with under optimal conditions
Application	IHC(P), ICC, ELISA

**Applications**

Somatotropin is a 25 kDa protein that can be expressed in the growth hormone receptor complex, extracellular space and extracellular region of cells. It is alternatively called Growth hormone 1, Pituitary growth hormone and Growth hormone. In humans, pituitary growth hormone reacts with growth hormone receptor, metal ion and prolactin receptor. Pituitary growth hormone has been closely linked to critical functions such as growth factor and hormone activity, and furthermore, is an integral part of organismal processes, for instance, animal organ development, response to estradiol and bone maturation. It is the subject of comprehensive examination because of the fact that it is involved with the growth hormone receptor signaling pathway via JAK-STAT, positive regulation of receptor signaling pathway via JAK-STAT and positive regulation of insulin-like growth factor receptor signaling pathway. Pituitary growth hormone promotes multicellular organism growth and MAP kinase activity. Mutations and abnormalities in pituitary growth hormone have been thought to be involved with a number of diseases, for example, isolated growth hormone deficiency 1B (IGHD1B), isolated growth hormone deficiency 1A (IGHD1A) and Kowarski syndrome (KWKS). Isolated growth hormone deficiency 1A- , an autosomal recessive inheritance disorder characterized by abnormality of the immune system, impaired growth-hormone response to insulin stimulation test and hypoglycemia, has in specific been of interest to scientists.