

**Purified Mouse Anti-human  $\alpha$ -Fetoprotein  
Antibody \*AFP-11, monoclonal\***

Catalog number: V103095

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	$\alpha$ -Fetoprotein
Clone	AFP-11

**Biological Properties**

---

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

**Applications**

---

$\alpha$ -fetoprotein (also known as  $\alpha$ -1-fetoprotein or  $\alpha$ -fetoglobulin) is a protein with a molecular weight of 69 kDa, found in the cytoplasm, endoplasmic reticulum lumen and extracellular space of cells. It is thought to be essential to organismal processes, in particular, ovulation from ovarian follicle, cellular protein metabolic process and progesterone metabolic process. Sequencing of  $\alpha$ -fetoprotein has shown it contains 3 conserved structural units: albumin 1, albumin 2 and albumin 3 domain.  $\alpha$ -fetoprotein binds with zinc ion and fatty acid. Mutations and abnormalities in  $\alpha$ -fetoprotein have been thought to be involved with a number of diseases, for instance,  $\alpha$ -fetoprotein, hereditary persistence (HPAFP) and  $\alpha$ -fetoprotein deficiency (AFPD).