

## Purified Mouse Anti-human Placental Alkaline Phosphatase Antibody \*H7E8, monoclonal\*

Catalog number: V1032025

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	Placental Alkaline Phosphatase
Clone	H7E8

### Biological Properties

---

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

### Applications

---

Alkaline phosphatase, placental type (alternatively called PLAP-1, Placental alkaline phosphatase 1 or Alkaline phosphatase Regan isozyme) is a transmembrane protein with a molecular weight of 58 kDa, located in the cell surface, plasma membrane and anchored component of membrane of cells. In humans, alkaline phosphatase Regan isozyme is an integral part of organismal processes, for example, dephosphorylation. Alkaline phosphatase Regan isozyme has been associated with important functions such as alkaline phosphatase activity. It recognizes GPI-anchor amidated aspartate, zinc ion and magnesium ion.