

**APC Mouse Anti-human/dog/pig TNAP
Antibody *W8B2B10, monoclonal***Catalog number: V1032340
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, dog, pig, sheep, monkey, goat
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	TNAP
Clone	W8B2B10
Conjugate	APC

Biological Properties

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

Applications

Alkaline phosphatase, tissue-nonspecific isozyme (AP-TNAP) is a 57 kDa protein that can be expressed in the anchored component of membrane, extracellular membrane-bounded organelle and plasma membrane of cells. In humans, AP-TNAP has been associated with key functions such as alkaline phosphatase and pyrophosphatase activity. AP-TNAP recognizes GPI-anchor amidated serine and metal ion. It plays an important role in organismal processes, namely, response to vitamin D, cellular response to organic cyclic compound and response to lipopolysaccharide. AP-TNAP is clinically significant because abnormalities in its function have been closely linked to diseases like hypophosphatasia (HOPS), hypophosphatasia childhood type (HOPSC) and hypophosphatasia infantile type (HOPSI). Hypophosphatasia infantile type, an autosomal recessive inherited disorder characterized by nephrocalcinosis, irritability and seizures, has in particular been of interest to scientists.