

---

# ATP12A Polyclonal Antibody

(Catalog # A72347)

---

## Background

The protein encoded by this gene belongs to the family of P-type cation transport ATPases. This gene encodes a catalytic subunit of the ouabain-sensitive H<sup>+</sup>/K<sup>+</sup> -ATPase that catalyzes the hydrolysis of ATP coupled with the exchange of H<sup>(+)</sup> and K<sup>(+)</sup> ions across the plasma membrane. It is also responsible for potassium absorption in various tissues. Two transcript variants encoding different isoforms have been found for this gene.

## Description

ATP12A Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

## Formulation

Buffer: PBS with 0.02% sodium azide, 50% glycerol, pH7.3.

## Specificity

Human

## Isotype

IgG

## Uniprot ID

P54707

## Purification

Affinity Purification

## Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 890-980 of human ATP12A (NP\_001172014.1).

## Storage

Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

## Alternative Names

ATP12A; ATP1AL1; HK; potassium-transporting ATPase alpha chain 2

## Application

WB; Recommended dilution: WB 1:500 - 1:2000

---