

## AQP1 Polyclonal Antibody

(Catalog A53228)

### Background

Forms a water-specific channel that provides the plasma membranes of red cells and kidney proximal tubules with high permeability to water, thereby permitting water to move in the direction of an osmotic gradient.

### Description

AQP1 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

### Formulation

Liquid. 0.03% Proclin 300, 50% Glycerol, 0.01M PBS, PH 7.4.

### Specificity

Human, Mouse, Rat

### Isotype

IgG

### Uniprot ID

P29972

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Human Aquaporin-1 protein (220-269AA)

### Storage

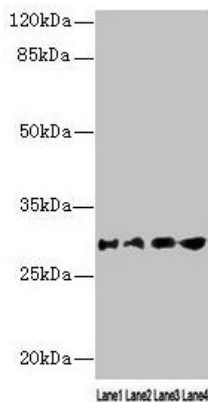
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

### Alternative Names

Aquaporin-CHIP, Urine water channel, Water channel protein for red blood cells and kidney proximal tubule, AQP1, CHIP28

### Application

ELISA, WB, IF, IP; Recommended dilution: WB:1:1000-1:5000, IF:1:50-1:200, IP:1:200-1:2000



#### Western blot

All lanes: AQP1 antibody at 16ug/ml

Lane 1: Rat heart tissue

Lane 2: Mouse skeletal muscle tissue

Lane 3: A549 whole cell lysate

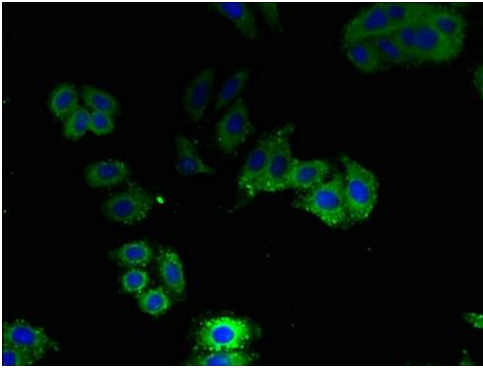
Lane 4: 293T whole cell lysate

Secondary

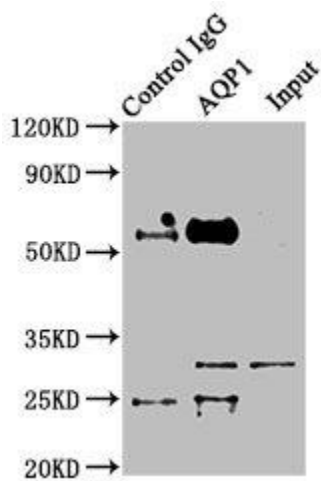
Goat polyclonal to rabbit IgG at 1/10000 dilution

Predicted band size: 29, 20, 24, 17 kDa

Observed band size: 29 kDa



Immunofluorescent analysis of HepG2 cells using AQP1 Antibody at dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)



Immunoprecipitating AQP1 in Mouse skeletal muscle tissue  
Lane 1: Rabbit control IgG instead of AQP1 Antibody in Mouse skeletal muscle tissue. For western blotting, a HRP-conjugated Protein G antibody was used as the secondary antibody (1/2000).  
Lane 2: AQP1 Antibody (8ug) + Mouse skeletal muscle tissue (500ug).  
Lane 3: Mouse skeletal muscle tissue (10ug).