

PVALB Polyclonal Antibody

(Catalog # A57009)

Background

In muscle, parvalbumin is thought to be involved in relaxation after contraction. It binds two calcium ions.

Description

PVALB Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

P20472

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Parvalbumin alpha protein (2-110AA)

Storage

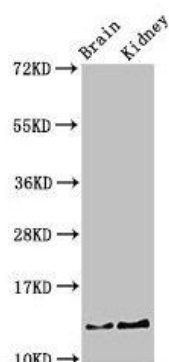
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

PVALB

Application

ELISA, WB, IHC, IF; Recommended dilution: WB 1:500 - 1:5000, IHC 1:20 - 1:200, IF 1:50 - 1:200



Western Blot

Positive WB detected in: Mouse brain tissue, Mouse kidney tissue

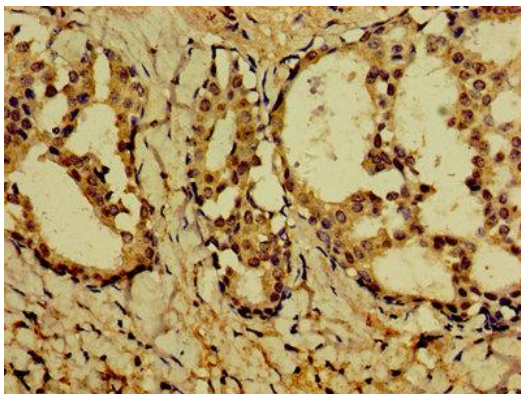
All lanes: PVALB Polyclonal Antibody at 3µg/ml

Secondary

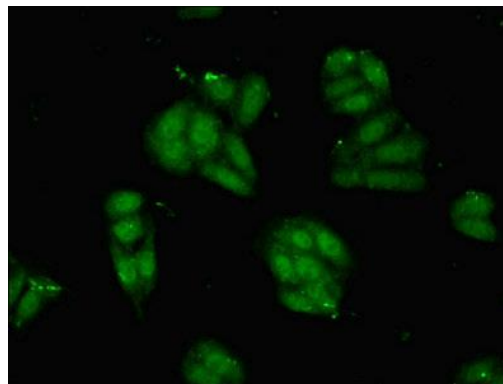
Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 13 kDa

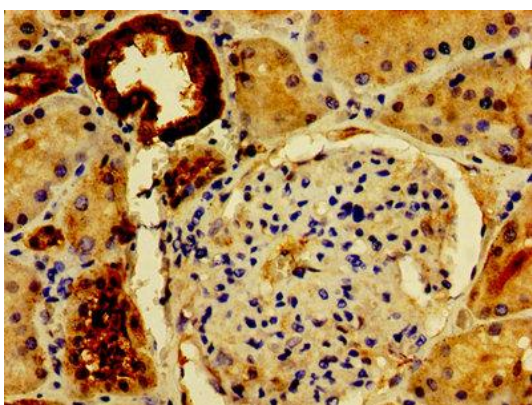
Observed band size: 13 kDa



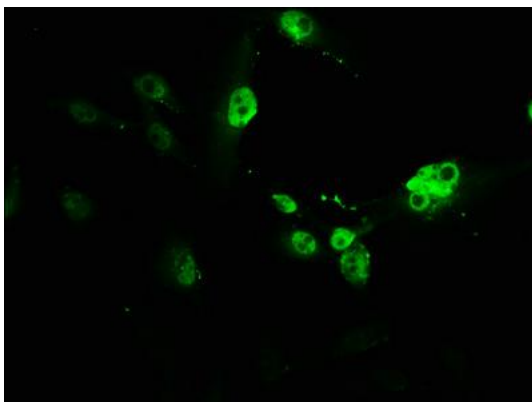
Immunohistochemistry of paraffin-embedded human breast cancer using PVALB Polyclonal Antibody at dilution of 1:100



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



Immunohistochemistry of paraffin-embedded human kidney tissue using PVALB Polyclonal Antibody at dilution of 1:100



Immunofluorescence staining of U251 cells with PVALB Polyclonal Antibody at 1:166, counter-stained with DAPI. The cells were fixed in 4% formaldehyde, permeabilized using 0.2% Triton X-100 and blocked in 10% normal Goat Serum. The cells were then incubated with the antibody overnight at 4°C. The secondary antibody was Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L).