

TRPV2 Polyclonal Antibody

(Catalog # A69355)

Background

is not activated by vanilloids and acidic pH.

Description

TRPV2 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

Q9Y5S1

Purification

Protein G purified

Immunogen

Recombinant Human Transient receptor potential cation channel subfamily V member 2 protein (1-117AA)

Storage

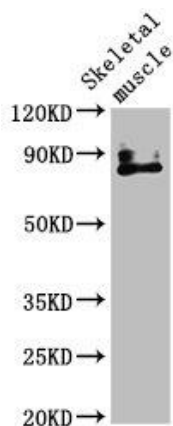
Shipped at 4°C. Upon receipt, store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Transient receptor potential cation channel subfamily V member 2, TrpV2, Osm-9-like TRP channel 2, OTRPC2, Vanilloid receptor-like protein 1, VRL-1, TRPV2, VRL

Application

ELISA, WB, IHC, IF; Recommended dilution: WB:1:500-1:5000, IHC:1:500-1:1000, IF:1:50-1:200



Western Blot

Positive WB detected in: Mouse skeletal muscle tissue

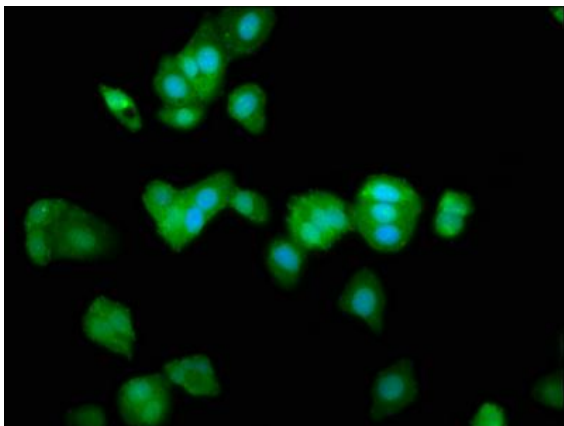
All lanes: TRPV2 antibody at 4.51 µg/ml

Secondary

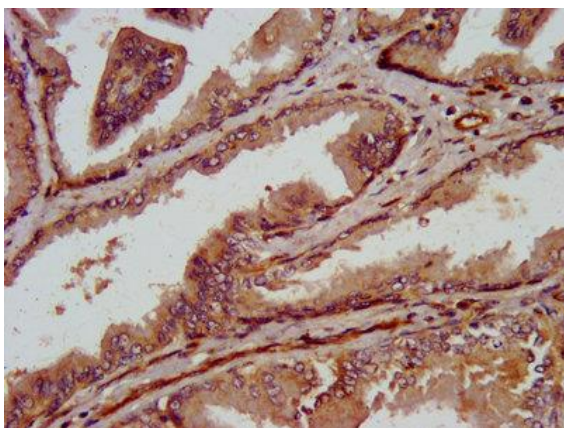
Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 86 kDa

Observed band size: 86 kDa



Immunofluorescence staining of HepG2 cells with TRPV2 Antibody at 1:266, 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).



IHC image of TRPV2 Antibody diluted at 1:800 and staining in paraffin-embedded human prostate tissue performed on a Leica Bond™ system. After dewaxing and hydration, antigen retrieval was mediated by high pressure in a citrate buffer (pH 6.0). Section was blocked with 10% normal goat serum 30min at RT. Then primary antibody (1% BSA) was incubated at 4°C overnight. The primary is detected by a biotinylated secondary antibody and visualized using an HRP conjugated SP system.