

# ASIC2 Polyclonal Antibody

(Catalog # A69851)

## Background

Cation channel with high affinity for sodium, which is gated by extracellular protons and inhibited by the diuretic amiloride. Also permeable for Li(+) and K(+). Generates a biphasic current with a fast inactivating and a slow sustained phase. Heteromeric channel assembly seems to modulate.

## Description

ASIC2 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

## Formulation

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

## Specificity

Human

## Isotype

IgG

## Uniprot ID

Q16515

## Purification

>95%, Protein G purified

## Immunogen

Recombinant Human Acid-sensing ion channel 2 protein (197-345AA)

## Storage

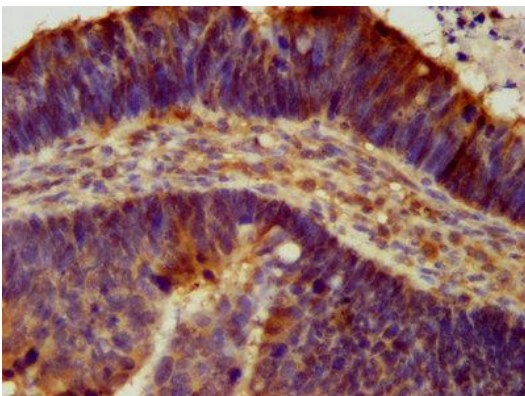
Shipped at 4°C. Upon receipt, store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

## Alternative Names

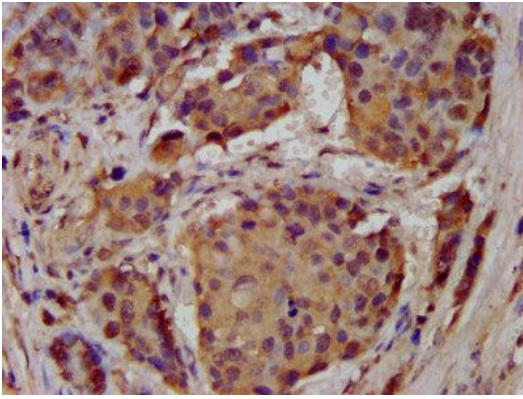
Acid-sensing ion channel 2, Amiloride-sensitive brain sodium channel, Amiloride-sensitive cation channel 1, neuronal, Amiloride-sensitive cation channel neuronal 1, Brain sodium channel 11, ASIC2, ACCN, ACCN1, BNAC1, MDEG

## Application

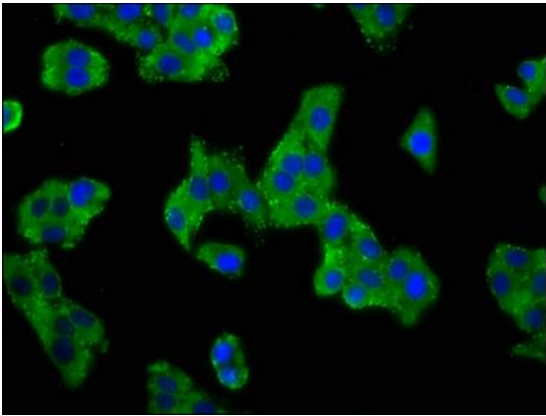
ELISA, IHC, IF; Recommended dilution: IHC: 1:200-1:500, IF: 1:50-1:200



IHC image of ASIC2 Polyclonal Antibody diluted at 1:300 and staining in paraffin-embedded human ovarian cancer 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.



IHC image of ASIC2 Polyclonal Antibody diluted at 1:300 and staining in paraffin-embedded human pancreatic cancer 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.



Immunofluorescence staining of HepG2 cells with ASIC2 Polyclonal Antibody at 1:100, 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).