
ASPH Polyclonal Antibody

(Catalog #A54361)

Background

Isoform 1: specifically hydroxylates an Asp or Asn residue in certain epidermal growth factor-like (EGF) domains of a number of proteins. Isoform 8: membrane-bound Ca^{2+} -sensing protein, which is a structural component of the ER-plasma membrane junctions. Isoform 8 regulates the activity of $\text{Ca}^{(+2)}$ released-activated $\text{Ca}^{(+2)}$ (CRAC) channels in T-cells.

Description

ASPH 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

Q12797

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Aspartyl/asparaginyl beta-hydroxylase protein (75-270AA)

Storage

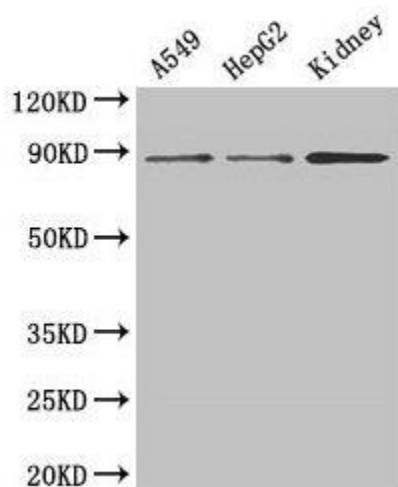
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

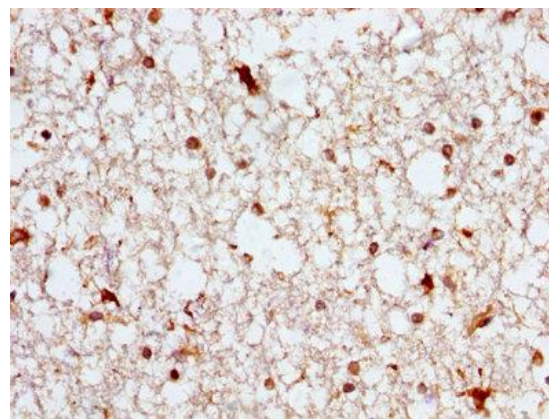
Aspartate beta-hydroxylase Peptide-aspartate beta-dioxygenase ASPH

Application

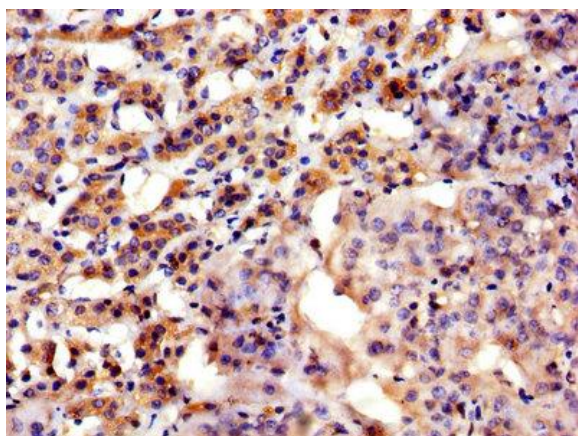
ELISA, WB, IHC, IF; Recommended dilution: WB:1:500-1:5000, IHC:1:200-1:500, IF:1:50-1:200



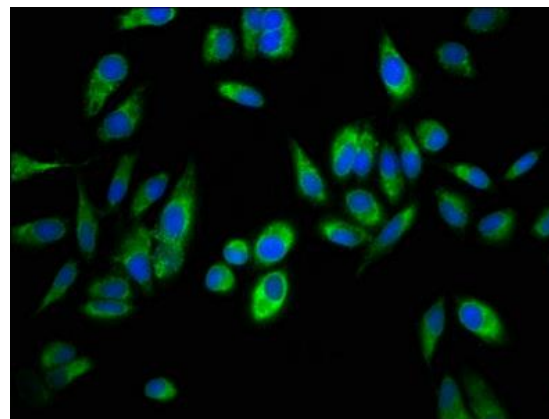
Western Blot
Positive WB detected in: A549 whole cell lysate, HepG2 whole cell lysate, Mouse kidney tissue
All lanes: ASPH Polyclonal Antibody at 3µg/ml
Secondary
Goat polyclonal to rabbit IgG at 1/50000 dilution
Predicted band size: 86, 35, 26, 24, 34, 84, 30, 22, 29, 33 kDa
Observed band size: 86 kDa



IHC image of ASPH Polyclonal Antibody diluted at 1:300 and staining in paraffin-embedded human brain 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.



IHC image of ASPH Polyclonal Antibody diluted at 1:300 and staining in paraffin-embedded human adrenal gland 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.



Immunofluorescence staining of HeLa cells with ASPH 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).