
SLC26A4 Polyclonal Antibody

(Catalog #A69707)

Background

Sodium-independent transporter of chloride and iodide.

Description

SLC26A4 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human

Isotype

IgG

Uniprot ID

O43511

Purification

Protein G purified

Immunogen

Recombinant Human Pendrin protein (592-655AA)

Storage

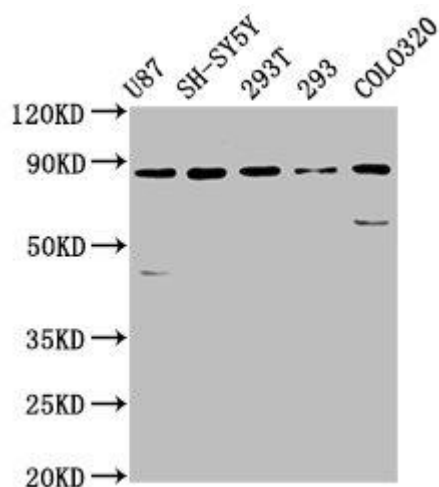
Shipped at 4°C. Upon receipt, store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Pendrin, Sodium-independent chloride/iodide transporter, Solute carrier family 26 member 4, SLC26A4, PDS

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: U87 whole cell lysate, SH-SY5Y whole cell lysate, 293T whole cell lysate, 293 whole cell lysate, COLO320 whole cell lysate

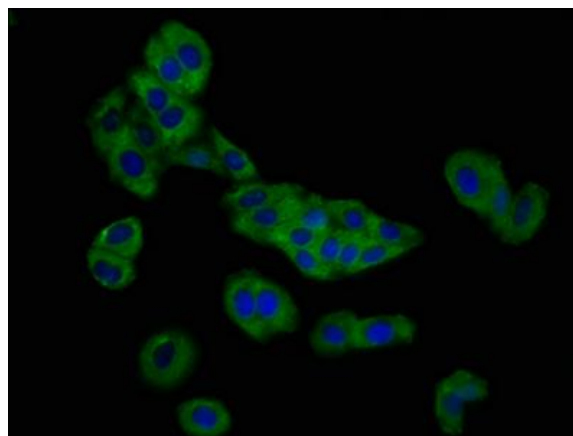
All lanes: SLC26A4 antibody at 5.7ug/ml

Secondary

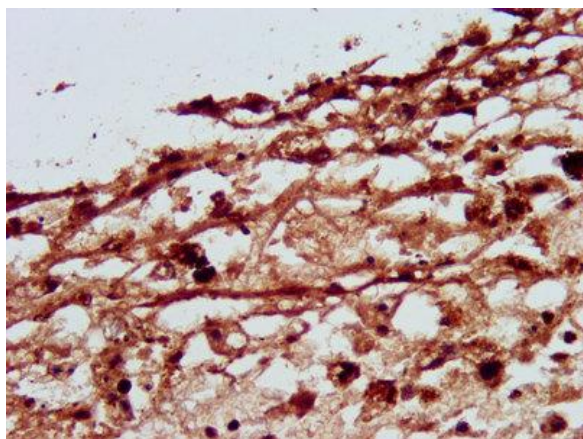
Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 86, 40 KDa

Observed band size: 86 KDa



Immunofluorescence staining of HepG2 cells with SLC26A4 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).



IHC image of SLC26A4 Polyclonal Antibody diluted at 1:500 and staining in paraffin-embedded human melanoma 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.