

SLC4A8 Polyclonal Antibody

(Catalog # A72220)

Background

The protein encoded by this gene is a membrane protein that functions to transport sodium and bicarbonate ions across the cell membrane. The encoded protein is important for pH regulation in neurons. The activity of this protein can be inhibited by 4,4'-Di-isothiocyanatostilbene-2,2'-disulfonic acid (DIDS). Several transcript variants encoding different isoforms have been found for this gene.

Description

SLC4A8 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Buffer: PBS with 0.02% sodium azide, 50% glycerol, pH7.3.

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

Q2Y0W8

Purification

Affinity Purification

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 1-90 of human SLC4A8 (NP_001035049.1).

Storage

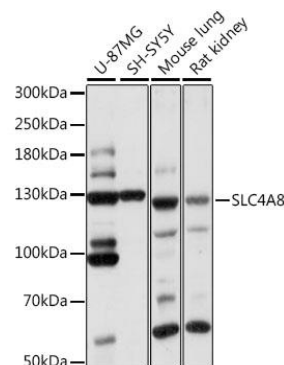
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

SLC4A8; NBC3; NDCBE; solute carrier family 4 member 8

Application

WB; Recommended dilution: WB 1:500 - 1:2000



Western blot analysis of extracts of various cell lines, using SLC4A8 Polyclonal Antibody at 1:1000 dilution. Secondary Antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution. Lysates/proteins: 25ug per lane. Blocking buffer: 3% nonfat dry milk in TBST. Exposure time: 1s.