
VAT1 Polyclonal Antibody

(Catalog #A73308)

Background

Synaptic vesicles are responsible for regulating the storage and release of neurotransmitters in the nerve terminal. The protein encoded by this gene is an abundant integral membrane protein of cholinergic synaptic vesicles and is thought to be involved in vesicular transport. It belongs to the quinone oxidoreductase subfamily of zinc-containing alcohol dehydrogenase proteins.

Description

VAT1 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Buffer: PBS with 0.01% thimerosal, 50% glycerol, pH7.3.

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

Q99536

Purification

Affinity Purified

Immunogen

A synthetic peptide corresponding to a sequence within amino acids 200-300 of human VAT1 (NP_006364.2).

Storage

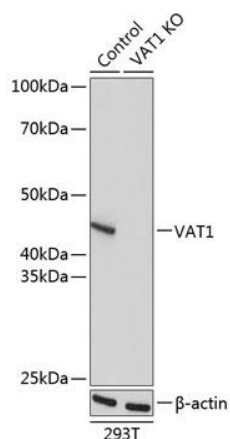
Shipped at 4°C. Store at -20°C. Avoid freeze / thaw cycles.

Alternative Names

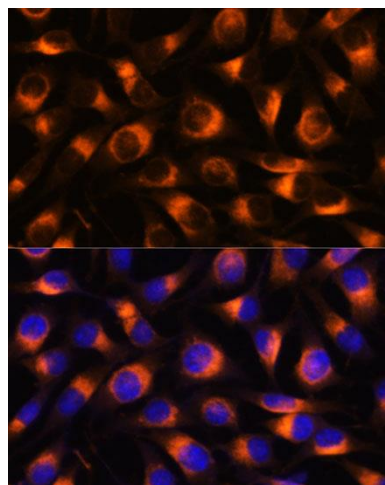
VAT1; VATI

Application

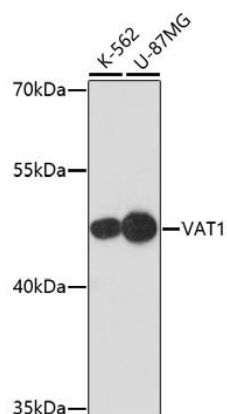
ELISA, WB, IF; Recommended dilution: WB 1:500 - 1:2000, IF 1:50 - 1:200



Western blot analysis of extracts from normal (control) and VAT1 knockout (KO) 293T cells, using VAT1 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.



Immunofluorescence analysis of L929 cells using VAT1 Polyclonal Antibody at dilution of 1:100. Blue: DAPI for nuclear staining.



Western blot analysis of extracts of various cell lines, using VAT1 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: 5min.