

MOLECULAR

**SIGNATURE®
VPS35
ELISA Kit**

Discovery through Partnership | Excellence through Quality

Colorimetric measurement of VPS35
Catalog No. SKT-141

Product Name

VPS35 ELISA Kit

Description

Colorimetric measurement of VPS35

Species Reactivity

Human, Mouse, Rat

Platform

Microplate

Sample Types

Cell Lysate, Tissue

Detection Method

Colorimetric

Assay Type

Sandwich ELISA (Enzyme-linked Immunosorbent Assay)

Sensitivity

50 pg/mL

Assay Range

78.125 - 5000 pg/mL

Incubation Time

3 hours

Number Of Samples

40 samples in duplicate

Other Resources

Kit Booklet Lot No. SV738133 , Kit Booklet Lot No. SV888883 , Kit Booklet Lot No. SV999550 , Kit Booklet Lot No. SV888770 , MSDS

Field Of Use

Not for use in humans. Not for use in diagnostics or therapeutics. For in vitro research use only.

Properties

Storage Temperature

4°C and -20°C

Shipping Temperature

Blue Ice

Product Type

ELISA Kits

Kit Overview

Component No.

Item

Quantity / Size

SKC-141A

Anti-VPS35 Immunoassay Plate

1 Plate

SKC-141B

5X VPS35 Extraction Reagent

1 bottle/10mL

SKC-141C

Recombinant VPS35 Standard

2 vials

SKC-141D

Standard and Sample Diluent (Red)

1 bottle/50mL

SKC-141E

10X Wash Buffer Concentrate

1 bottle/100mL

SKC-141F

Anti-VPS35 Biotinylated Antibody Concentrate

1 vial/150uL

SKC-141G

Anti-VPS35 Biotinylated Antibody Diluent (Green)

1 bottle/13mL

SKC-141H

100X Streptavidin:HRP Concentrate

1 vial/150uL

SKC-141I

Streptavidin: HRP Diluent (Purple)

1 bottle/13mL

SKC-141J

TMB Substrate

1 bottle/13mL

SKC-141K

Stop Solution

1 bottle/13mL

Biological Description

Alternative Names

FLJ10752, MEM3, PARK17, TCCCTA00141, VPS35 retromer complex component, Vacuolar protein sorting-associated protein 35, maternal-embryonic 3, vesicle protein sortin 35

Research Areas

Alzheimer's Disease, Cell Signaling, Golgi Proteins, Membrane Trafficking Proteins, Neurodegeneration, Neuroscience, Parkinson's Disease, Protein Trafficking

Scientific Background

Vacuolar Protein Sorting 35 (VPS35) is a core component of the retromer complex, which mediates the recycling of membrane proteins from endosomes to the Golgi apparatus. VPS35 plays a critical role in maintaining neuronal protein homeostasis, synaptic function, and mitochondrial integrity.

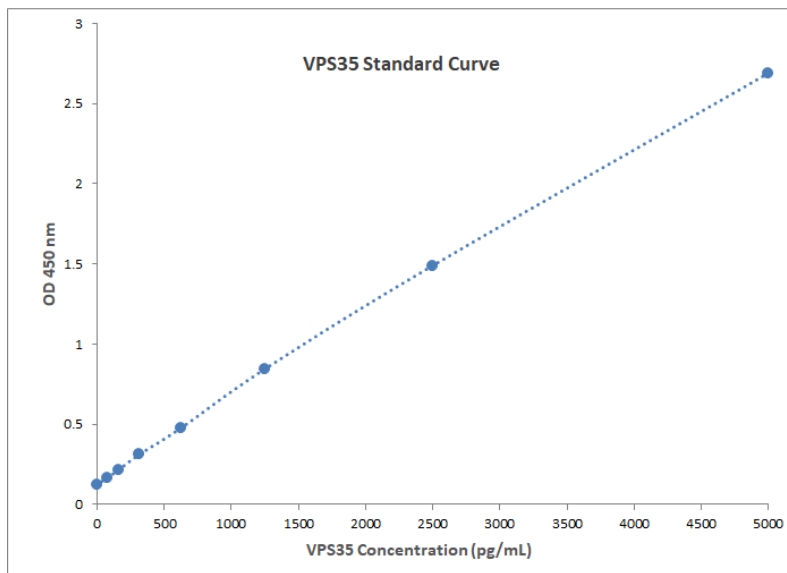
Mutations in VPS35, particularly the D620N variant, have been directly linked to familial Parkinson's disease. These mutations impair retromer function, leading to defective trafficking of key neuronal receptors and enzymes, increased protein aggregation, and mitochondrial fragmentation.

In neurodegeneration research, VPS35 is a focal point for understanding how intracellular transport defects contribute to dopaminergic neuron loss. Targeting retromer dysfunction offers a promising therapeutic strategy for modifying disease progression in Parkinson's and related disorders.

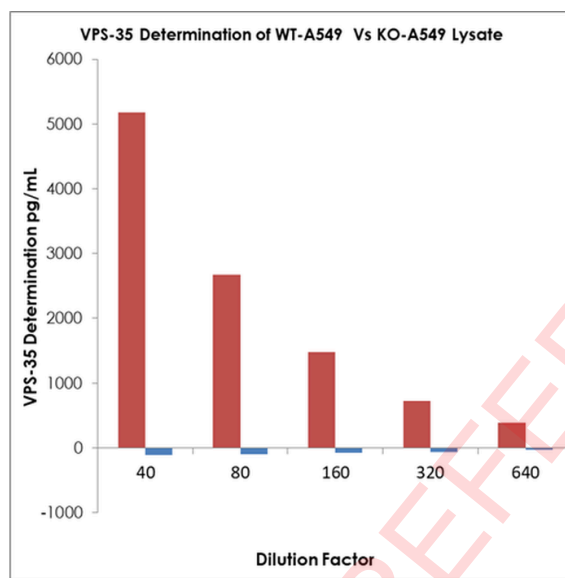
References

1. Vilarino-Guell, C. et al. (2011) Am J Hum Genet 89:162–167
 2. Zimprich, A. et al. (2011) Am J Hum Genet 89:168–175
 3. Rahman, A.A., Morrison, B.E. (2019) Neurosci 401:1-10.
-

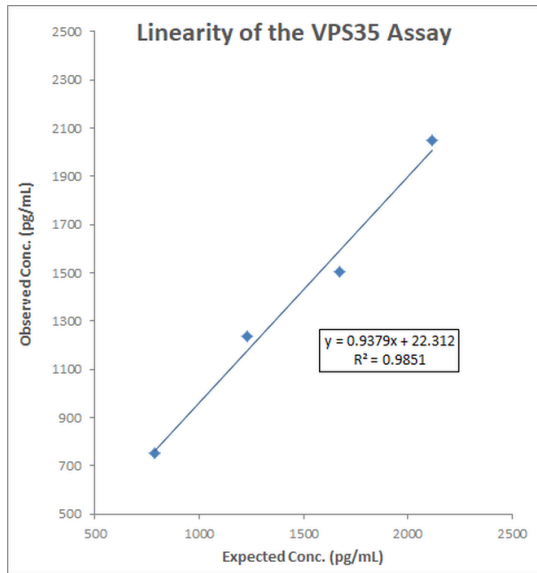
Product Images



Typical Standard Curve for the VPS35 Kit (Enzyme-Linked Immunosorbent Assay) MOLECULAR SIGNATURE® – SKT-141. Assay Type: Competitive ELISA. Detection Method: Colorimetric Assay.



VPS35 Determination of WT-A549 vs KN-549 Lysates Graph for the VPS35 Kit (Enzyme-Linked Immunosorbent Assay) MOLECULAR SIGNATURE® – SKT-141. Assay Type: Competitive ELISA. Detection Method: Colorimetric Assay.



Linearity was determined by using low (rat kidney) and high (mouse kidney) natural samples, diluted 1/100 and mixed in various proportions.

Reviews

There are no reviews yet.

Product Citations
