

STK25 Polyclonal Antibody

(Catalog # A71397)

Background

This gene encodes a member of the germinal centre kinase III (GCK III) subfamily of the sterile 20 superfamily of kinases. The encoded enzyme plays a role in serine-threonine liver kinase B1 (LKB1) signaling pathway to regulate neuronal polarization and morphology of the Golgi apparatus. The protein is translocated from the Golgi apparatus to the nucleus in response to chemical anoxia and plays a role in regulation of cell death. A pseudogene associated with this gene is located on chromosome 18. Multiple alternatively spliced transcript variants have been observed for this gene.

Description

STK25 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

O00506

Purification

Affinity Purification

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 297-426 of human STK25 (NP_006365.2).

Storage

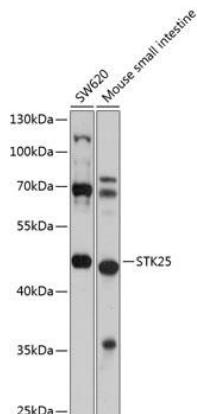
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

STK25; SOK1; YSK1; serine/threonine kinase 25

Application

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



Western blot analysis of extracts of various cell lines, using STK25 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: 15s.