

Anti-TTBK1

Rabbit Polyclonal Antibody

Catalog # T18-63R

Lot # O2121-19

Cited Applications

WB, ELISA, IF, IHC

Ideal working dilutions for each application should be empirically determined by the investigator.

Specificity

Recognizes the TTBK1 protein

Cross Reactivity

Human and Mouse

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

TTBK1 antibody was raised against a 16 amino acid synthetic peptide near the carboxy terminus of human TTBK1

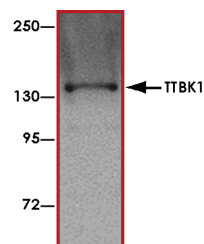
Formulation

PBS + 0.02% sodium azide

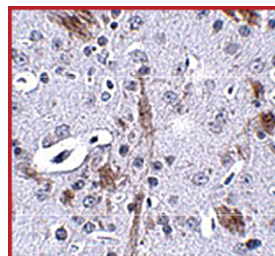
Stability

1yr at -20°C from date of shipment

Sample Data



Western blot analysis of TTBK1 in Jurkat lysate with TTBK1 antibody at 1 µg/ml.



Immunohistochemistry of TTBK1 in mouse brain tissue with TTBK1 antibody at 2.5 µg/ml.

Anti-TTBK1

Rabbit Polyclonal Antibody

Catalog Number

T18-63R

Specific Lot Number

O2121-19

Purification

Affinity chromatography

Stability

1yr at -20°C from date of shipment

Storage & Shipping

Store product at -20°C. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on ice packs.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.