

## Anti-Stathmin

Rabbit Polyclonal Antibody

**Catalog # S71-363R**

Lot # O2121-7

### Cited Applications

WB, ELISA, IF, IHC

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

### Specificity

Recognizes the Stathmin protein

### Cross Reactivity

Human and Mouse

### Host/Isotype/Clone#

Rabbit, IgG

### Immunogen

Stathmin antibody was raised against a 20 amino acid synthetic peptide from near the carboxy terminus of human Stathmin

### Formulation

PBS + 0.02% sodium azide

### Stability

1yr at -20°C from date of shipment

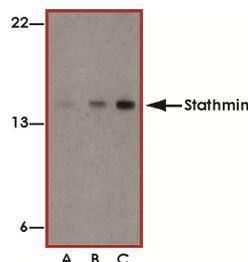
### Scientific Background

Stathmin is a ubiquitous, highly conserved protein present in the cytoplasm in a variety of unphosphorylated and phosphorylated forms. Stathmin functions as an intracellular relay integrating regulatory signals of the cellular environment (1). Stathmin is involved in the regulation of the microtubule filament system. Stathmin binds tubulin in a ternary complex, with 2 tubulins for every stathmin and this complex interferes with microtubule dynamics in vitro and in vivo. Overexpression of several stathmin mutants suggest that stathmin induces depolymerization of interphase and mitotic microtubules in its unphosphorylated state, but is inactivated by phosphorylation in mitosis (2).

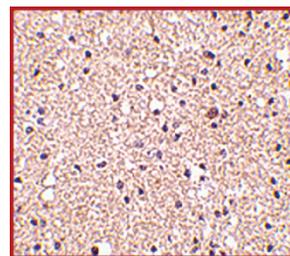
### References

1. Sobel, A. Stathmin: a relay phosphoprotein for multiple signal transduction? Trends Biochem. Sci. 16: 301-305, 1991.
2. Gavet, O. et al: The stathmin phosphoprotein family: intracellular localization and effects on the microtubule network. J. Cell Sci. 111: 3333-3346, 1999.

### Sample Data



Western blot analysis of Stathmin in EL4 cell lysate with Stathmin antibody at (A) 0.5, (B) 1 and (C) 2 ug/ml.



Immunohistochemistry of Stathmin in human brain tissue with Stathmin antibody at 2.5 ug/mL.

## Anti-Stathmin

Rabbit Polyclonal Antibody

Catalog Number

S71-363R

Specific Lot Number

O2121-7

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](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

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