

## Anti-ELK1

Rabbit Polyclonal Antibody

### Catalog # E64-63R

Lot # M191-4

### Cited Applications

- Western blot (1:1000)

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

### Specificity

Recognizes the ELK1 protein

### Cross Reactivity

- Western blot human ELK1

ELK1 from other species may also be detectable

### Host

Rabbit

### Immunogen

Synthetic peptide corresponding to amino acids 387-391 of human ELK1

### Formulation

PBS, pH7.4, 0.02%NaN<sub>3</sub>, 50% glycerol

### Stability

Store at 4°C (add 0.1% NaN<sub>3</sub>) for several months, and at -20°C for longer periods. 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.

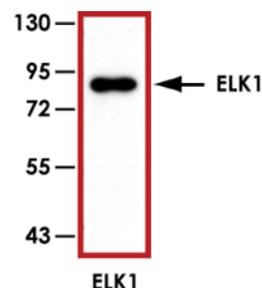
### Scientific Background

ELK1 is a member of the ETS oncogene family of transcription factors (1). ELK1 protein interacts with promoter of the c-fos proto-oncogene and has been shown to be involved in the Ras-Raf-MAPK signaling cascade (1). Immature phosphorylation of transcription factor ELK-1 is implicated in premature aging syndrome and insulin resistance. Expression studies have revealed that elevated expression of ELK1-like protein in human esophageal can result in squamous cell carcinoma (2).

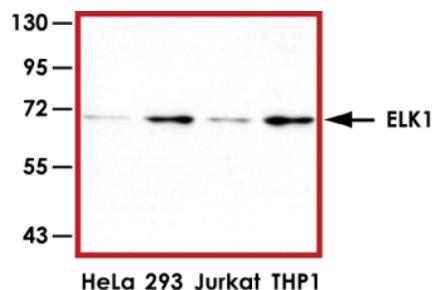
### References

1. Chen, A.G. et al: Overexpression of Ets-like protein 1 in human esophageal squamous cell carcinoma. *World J Gastroenterol.* 2006, 28;12(48):7859-63.
2. Knebel, B. et al: Reduced phosphorylation of transcription factor Elk-1 in cultured fibroblasts of a patient with premature aging syndrome and insulin resistance. *Exp Clin Endocrinol Diabetes.* 2005;113(2):94-101.

### Sample Data



Representative Western blot with Anti-ELK1 (1:1000) using 15ng of GST-tagged human recombinant ELK1.



Representative Western blot with Anti-ELK1 (1:1000) using 20µg of HeLa, 293, Jurkat and THP1 cell lysates.

## Anti-ELK1

Rabbit Polyclonal Antibody

Catalog Number

E64-63R

Specific Lot Number

M191-4

Purification

Affinity chromatography

Concentration

1.0 µg/µL

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.**