

## Elk-1 antibody (pAb)

**Catalog Nos:** 61301, 61302

**RRID:** AB\_2793587

**Isotype:** IgG

**Application(s):** IP, WB

**Reactivity:** Human

**Volumes:** 100 µl, 10 µl

**Purification:** Affinity Purified

**Host:** Rabbit

**Molecular Weight:** 68 kDa

**Background:** Elk-1 (ETS-like gene 1) is a member of the ETS family of transcription factors and stimulates transcription. It binds to purine-rich DNA sequences, and can form a ternary complex with the serum response factor and the ETS and SRF motifs of the fos serum response element.

**Immunogen:** This Elk-1 antibody was raised against a peptide within the C-terminal region of human Elk-1.

**Buffer:** Purified IgG in PBS with 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

### Application Notes:

Applications Validated by Active Motif:

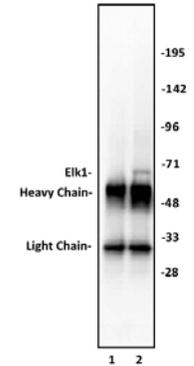
IP: 10 µl per IP

WB: 1:500 - 1:1,000 dilution

The addition of 0.05% Tween 20 in the blocking buffer and primary antibody incubation buffer is recommended to aid in detection by Western blot. Individual optimization may be required.

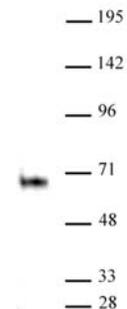
**Storage and Guarantee:** Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



### Elk-1 antibody (pAb) tested by Immunoprecipitation.

10 µl of Elk-1 antibody was used to immunoprecipitate Elk-1 from 400 µg of HeLa whole cell extract (lane 2). 10 µl of rabbit IgG was used as a negative control (lane 1). The immunoprecipitated protein was detected by Western blotting using the Elk-1 antibody at a dilution of 1:500.



### Elk-1 antibody (pAb) tested by Western blot.

Whole cell extract of HeLa cells (40 µg) probed with Elk-1 antibody (1:500).