

LexA DNA-binding Domain antibody (pAb)

Catalog Nos: 39184, 39185**RRID:** AB_2793174**Isotype:** Serum**Application(s):** ChIP, WB**Reactivity:** Budding Yeast, Not Species Specific**Volumes:** 100 µl, 10 µl**Purification:** None**Host:** Rabbit

Background: LexA is a prokaryotic protein that normally functions to repress the expression of DNA-damage response genes. The LexA DNA binding domain is frequently used as an epitope tag or fused to the "bait" protein as part of the yeast two-hybrid protein interaction system.

Immunogen: This LexA DNA-binding Domain antibody was raised against recombinant protein corresponding to full-length *E. coli* LexA.

Buffer: Rabbit serum containing 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

Applications Validated by Active Motif:

IP: 1 µl/ 200 µg lysate

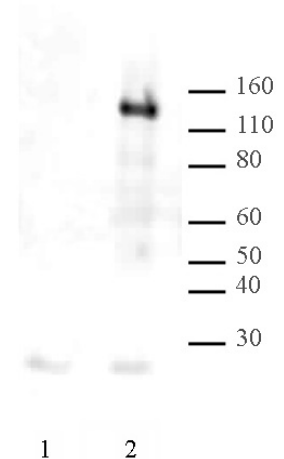
WB*: 1:25,000 - 1:100,000 dilution

Validation for Western blot was confirmed using a LexA-Rb fusion and a LexA-PP2A, β subunit fusion expressed in budding yeast. Individual optimization may be required.

*Note: many chromatin-bound proteins are not soluble in a low salt nuclear extract and fractionate to the pellet. Therefore, we recommend a High Salt / Sonication Protocol when preparing nuclear extracts for Western blot.

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.

**LexA DNA-binding Domain pAb tested by Western blot.**

20 µg of crude yeast protein extracts probed with LexA DNA-binding Domain pAb (1:100,000 dilution).

Lane 1: Extract from wild-type yeast.

Lane 2: Extract from yeast expressing a LexA-Retinoblastoma (Rb) fusion protein.