

SALL4 antibody (pAb)

Catalog Nos: 39957, 39958

RRID: AB_2793408

Isotype: IgG

Application(s): IP, WB

Reactivity: Mouse

Volumes: 100 μ l, 10 μ l

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 142 kDa

Background: SALL4 (Sal-like protein 4) is a zinc-finger transcription factor required for the maintenance of embryonic stem cell pluripotency. SALL4 activates the expression of Oct4 and reduction in the levels of SALL4 eliminates the ability of ES cells to self-renew and causes differentiation. SALL4 mutations are associated with the developmental disorders IVIC syndrome and Duane-radial ray syndrome.

Immunogen: This SALL4 antibody was raised against a synthetic peptide derived from the N-terminus of human SALL4.

Buffer: Purified IgG in 70 mM Tris (pH 8), 105 mM NaCl, 31 mM glycine, 0.07 mM EDTA, 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

Applications Validated by Active Motif:

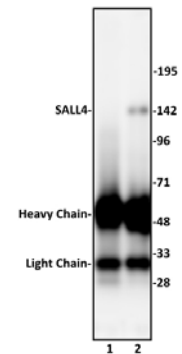
IP: 10 μ l per IP

WB*: 1:500 - 1:2,000 dilution

*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.

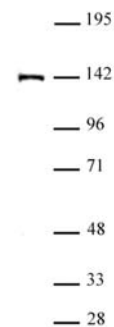
Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product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

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



SALL4 antibody (pAb) tested by Immunoprecipitation.

10 μ l of SALL4 antibody was used to immunoprecipitate SALL4 from 500 μ g of F9 nuclear cell extract (lane 2). 10 μ l of rabbit IgG was used as a negative control (lane 1). The immunoprecipitated protein was detected by Western blotting using the SALL4 antibody at a dilution of 1:500.



SALL4 antibody (pAb) tested by Western blot.

Nuclear extract of P19 cells (20 μ g) probed with SALL4 antibody (1:500).