

GLI1 antibody (pAb)

Catalog Nos: 61215, 61216

RRID: AB_2614953

Isotype: IgG

Application(s): WB

Reactivity: Human

Volumes: 100 μ l, 10 μ l

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 120 kDa

Background: GLI1 (Glioma-associated oncogene 1) is a zinc finger transcription factor and oncogene. GLI1 is part of the hedgehog signaling pathway, but is also activated in response to TGF- β signaling as well. GLI1 activates transcription of a number of genes that promote cell cycle progression. GLI1 is overexpressed as a result of constitutive activation of hedgehog signaling, which plays a role in the initiation of cancer. GLI1 also promotes self-renewal of neural stem cells.

Immunogen: This GLI1 antibody was raised against a peptide within the C-terminal region of human GLI1.

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

Application Notes:

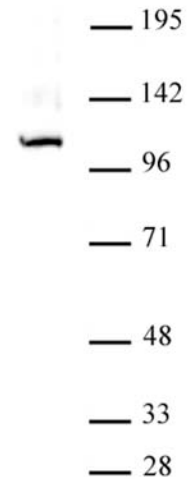
Applications Validated by Active Motif:

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.



GLI1 antibody (pAb) tested by Western blot.

Nuclear extract of Raji cells (20 μ g) probed with GLI1 antibody (pAb) at a dilution of 1:1,000.