

STAT3 antibody (pAb)

Catalog Nos: 61645, 61646

RRID: AB_2793718

Isotype: IgG

Application(s): WB

Reactivity: Human

Volumes: 100 μ l, 10 μ l

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 96 kDa

Background: **STATs** (signal transducers and activators of transcription) are a family of 7 transcription factors that form part of the JAK-STAT signaling cascade. This cascade is the basis of the signal transduction mechanism for many cytokine receptors. Once activated by phosphorylation by JAKs, STATs translocate to the nucleus. Accumulation of STATs in the nucleus is both rapid and tightly controlled. A number of factors regulate the JAK-STAT pathway including STAT dephosphorylation by phosphatases, altered nuclear import-export dynamics of STAT, and STAT gene activation antagonists such as SOCS (suppressors of cytokine signaling) and PIAS (Protein Inhibitors of Activated STATs).

Immunogen: This antibody was raised against a peptide within the N-terminal region of human STAT3.

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

Application Notes:

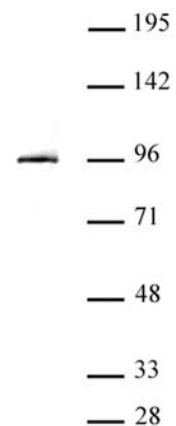
Applications Validated by Active Motif:

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

*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.



STAT3 antibody (pAb) tested by Western blot.

Detection of STAT3 by Western blot analysis. Nuclear extract of HepG2 cells (30 μ g) probed with STAT3 antibody at a dilution of 1:2,000.