

STAT3 phospho Ser727 antibody (pAb)

Catalog Nos: 39613, 39614

RRID: AB_2793276

Isotype: IgG

Application(s): DB, WB

Reactivity: Human

Volumes: 200 µl, 10 µl

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 92 kDa

Background: STAT3 (signal transducer and activator of transcription 3) is a member of the STAT family of transcription factors. STATs undergo tyrosine and serine phosphorylation in response to growth factor or cytokine signaling. This phosphorylation results in dimerization and translocation of STAT proteins to the nucleus. The STAT3 protein is expressed in most cell types and is activated by multiple cytokines and growth factors including: IFN α , IL-10, IL-6, IL-11, IL-12, IL-2, EGF, among others.

Immunogen: This STAT3 phospho Ser727 antibody was raised against a peptide containing phospho Ser727 of human STAT3.

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:

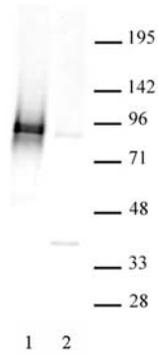
Applications Validated by Active Motif:

WB: 1:1,000 - 1:5,000 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.



STAT3 phospho Ser727 pAb tested by Western blot.

Hep G2 nuclear extract (15 µg per lane) was probed with STAT3 phospho Ser727 pAb (1:1,000).

Lane 1: cells stimulated with IL-6 (100 ng/ml for 15 minutes at 37°C).

Lane 2: No treatment.

STAT3 phospho Ser727 pAb tested by dot blot analysis.

Dot blot analysis was used to confirm the specificity of STAT3 phospho Ser727 pAb for STAT2 phospho Ser727. Phosphorylated peptides corresponding to the immunogen and related peptides were spotted onto PVDF and probed with the antibody at 1:30,000. The amount of peptide (picomoles) spotted is indicated next to each row.



Lane 1: Unmodified Ser727 STAT1 peptide.

Lane 2: Phospho Ser727 STAT1 peptide.

Lane 3: Unmodified Tyr689 STAT2 peptide.

Lane 4: Phospho Tyr689 STAT2 peptide.

Lane 5: Unmodified Ser727 STAT3 peptide.

Lane 6: Phospho Ser727 STAT3 peptide.

Lane 7: Unmodified Tyr705 STAT3 peptide.