

STAT5A/B phospho Tyr694/Tyr699 antibody (pAb)

Catalog Nos: 39617, 39618

RRID: AB_2793277

Isotype: IgG

Application(s): DB, WB

Reactivity: Human

Volumes: 200 µl, 10 µl

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 92 kDa

Background: STAT5A/B (signal transducer and activator of transcription 5) is a member of the STAT family of transcription factors. STATs undergo serine and tyrosine phosphorylation in response to growth factor or cytokine signaling. This phosphorylation results in dimerization and translocation of STAT proteins to the nucleus. The STAT5 protein is expressed in most cell types and is activated by multiple cytokines and growth factors including: IFNA1, IL-10, IL-6, IL-11, IL-12, IL-2, EGF, among others. Other members of the STAT family include STAT1 α , STAT2, STAT3, STAT4, STAT5A, STAT5B and STAT6.

Immunogen: This STAT5A/B phospho Tyr694/Tyr699 antibody was raised against a peptide containing phospho Tyr694 of human STAT5A. Human STAT5B shares the identical sequence surrounding Tyr699.

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.

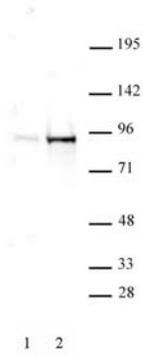
Application Notes:

Applications Validated by Active Motif:

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

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.



STAT5A/B phospho Tyr694/Tyr699 rabbit pAb tested by Western blot.

Nb2 nuclear extract (20 µg per lane) was probed with STAT5A/B phospho Tyr694/Tyr699 rabbit pAb (1:2,000).

Lane 1: No treatment.

Lane 2: Cells treated with prolactin (25 ng/ml for 20 minutes at 37°C).

STAT5A/B phospho Tyr694/Tyr699 rabbit pAb tested by dot blot analysis.

Dot blot analysis was used to confirm the specificity of STAT5A/B phospho Tyr694/Tyr699 rabbit pAb for STAT5A/B phospho Tyr694/699. 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. Lane 8: Phospho Tyr705 STAT3 peptide. Lane 9: Unmodified Ser726 STAT5A/Ser731 STAT5B peptide. Lane 10: Phospho Ser726 STAT5A/Ser731 STAT5B peptide. Lane 11: Unmodified Tyr694 STAT5A/Tyr699 STAT5B peptide. Lane 12: Phospho Tyr694 STAT5A/Tyr699 STAT5B peptide.