

STAT2 phospho Tyr689 antibody (pAb)

Catalog Nos: 39611, 39612

RRID: AB_2793275

Isotype: IgG

Application(s): DB, WB

Reactivity: Human

Volumes: 200 µl, 10 µl

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 110 kDa

Background: STAT2 (signal transducer and activator of transcription 2) 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. STAT2 has been shown to be activated by IFN α and IFN β , and to associate with the transcription factors p48 and p300.

Immunogen: This STAT2 phospho Tyr689 antibody was raised against a peptide containing phospho Tyr689 of human STAT2.

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.

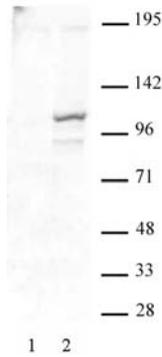
Application Notes:

Applications Validated by Active Motif:

WB: 1:500 - 1:1,000 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.



STAT2 phospho Tyr689 pAb tested by Western blot.

HeLa whole-cell extract (20 µg per lane) was probed with STAT2 phospho Tyr689 pAb (1:500).

Lane 1: No treatment.

Lane 2: Cells treated with IFNα (2500 units/ml for 30 minutes).

STAT2 phospho Tyr689 pAb tested by dot blot analysis.

Dot blot analysis was used to confirm the specificity of STAT2 phospho Tyr689 pAb for STAT2 phospho Tyr689. Phosphorylated peptides corresponding to the immunogen and related peptides were spotted onto PVDF and probed with the antibody at 1:10,000. 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.