

STAT4 antibody (pAb)

Catalog Nos: 61507, 61508

RRID: AB_2793663

Isotype: IgG

Application(s): WB

Reactivity: Human, Mouse

Volumes: 100 μ l, 10 μ l

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 86 kDa

Background: STAT4 (Signal Transducer And Activator Of Transcription 4) is involved in gene activation in response to IL-12, and is required for the development of the Th1 cell lineage. Both IL-12 and IFN α are capable of activating STAT4. Variants of the STAT4 gene are frequently associated with numerous autoimmune disorders.

Immunogen: This antibody was raised against a peptide within the C-terminal region of human STAT4.

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

Application Notes:

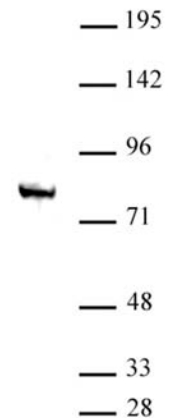
Applications Validated by Active Motif:

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

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.



STAT4 antibody (pAb) tested by Western blot.

Nuclear extract of mouse testis (30 μ g) probed with STAT4 antibody (pAb) at a dilution of 1:2,000.