

IRF-5 antibody (pAb)

Catalog Nos: 61269, 61270

RRID: AB_2615035

Isotype: IgG

Application(s): ChIP, ChIP-Seq, WB

Reactivity: Human

Volumes: 100 µl, 10 µl

Purification: Affinity Purified

Host: Rabbit

Molecular Weight: 60 kDa

Background: IRF-5 (interferon regulatory factor 5) is a transcription factor involved in the induction of interferons IFNA and INFB and inflammatory cytokines upon virus infection. Activated by TLR7 or TLR8 signaling.

Immunogen: This antibody was raised against a peptide within the N-terminal region of human IRF-5.

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:

ChIP: 10 µl per ChIP

ChIP-Seq & ChIP-chip: 10 µl each

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

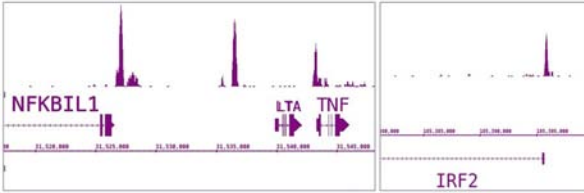
The addition of 0.05% Tween 20 in the blocking buffer and primary antibody incubation buffer is recommended to aid in detection by Western blot. Individual optimization may be required.

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.

IRF-5 antibody (pAb) tested by ChIP-Seq.

(ChIP) was performed using the ChIP-IT[®] High Sensitivity Kit (Cat. No. 53040) and 25 µg of chromatin from primary human macrophages + compound. ChIP DNA was sequenced on the Illumina HiSeq and 8 million sequence tags were mapped to identify IRF-5 binding sites across the genome. Both images show the expected enrichment.



IRF-5 antibody (pAb) tested by Western blot

Nuclear extract of THP-1 cells (20 µg) probed with IRF-5 antibody (1:500).

