

IRF-3 antibody (pAb)

Catalog Nos: 39371, 39372

RRID: AB_2793232

Isotype: Serum

Application(s): ChIP, WB

Reactivity: Human

Volumes: 200 μ l, 10 μ l

Purification: None

Host: Rabbit

Molecular Weight: 50 kDa

Background: IRF-3 (Interferon regulatory factor 3 or IRF3) is a member of the IRF family of transcription factors that include IRF-1, IRF-6 and IRF-7. IRF-3 that has been phosphorylated at serine/threonine forms a complex with CREBBP and translocates to the nucleus to activate transcription of interferon-induced genes.

Immunogen: This IRF-3 antibody was raised against a peptide corresponding to residues 306-312 of human IRF-3.

Buffer: Rabbit serum containing 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

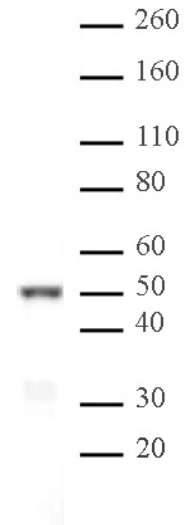
Application Notes:

Applications Validated by Active Motif:

WB: 1:1,000 - 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.



IRF-3 pAb tested by Western blot.

Detection of IRF-3 by Western blot analysis. The presence of IRF-3 in nuclear extracts derived from U-937 (IFN γ) cells is confirmed using IRF-3 rabbit polyclonal antibody at a 1:1,000 dilution.