

## IRAK-2 antibody (pAb)

**Catalog No:** 40912

**RRID:** AB\_2793434

**Application(s):** WB

**Reactivity:** Human

**Quantity:** 100 µg

**Purification:** Affinity Purified

**Host:** Rabbit

**Isotype:** IgG

**Concentration:** 0.5 µg/µl

**Molecular Weight:** 65 kDa

**Background:** IRAK-2 – Interleukin-1 Receptor-Associated Kinase2 is one of two serine/threonine kinases that become associated with the interleukin-1 receptor (IL1R) upon stimulation to upregulate NFκB (NFκB p50 & NFκB p65) expression.

**Immunogen:** This IRAK-2 antibody was raised against a peptide corresponding to amino acid residues 546-564 of human IRAK-2.

**Buffer:** PBS containing 0.02% sodium azide. Sodium azide is highly toxic.

**Application Notes:**

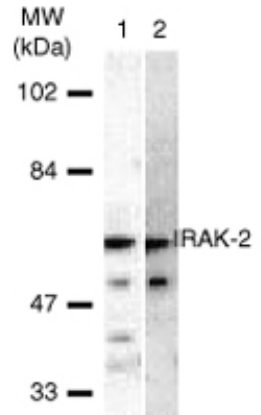
Applications Validated by Active Motif:

WB: 1 - 2 µg/ml dilution

For optimal results, primary antibody incubations should be performed at room temperature. The addition of 0.1% Tween 20 to all blocking solutions may also reduce background. Individual optimization may be required.

**Storage and Guarantee:** Some products may be shipped at room temperature. This will not affect their stability or performance. Store at 4°C for short term. 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.



**IRAK-2 pAb tested by Western blot.**

Detection of IRAK-2 by Western blot analysis. The analysis was performed using 10 µg/lane of HeLa (lane 1) and K-562 (lane 2) nuclear extracts and IRAK-2 pAb at a 2 µg/ml dilution.