

## EGF Receptor phospho Tyr992 antibody (pAb)

**Catalog No:** 39094

**RRID:** AB\_2793153

**Isotype:** IgG

**Application(s):** WB

**Reactivity:** Human

**Volume:** 200 µl

**Purification:** Affinity Purified

**Host:** Rabbit

**Concentration:** 0.05 µg/µl

**Molecular Weight:** 170 kDa

**Background:** EGFR – Epidermal Growth Factor Receptor is a member of the superfamily of transmembrane receptor tyrosine kinases and is widely expressed. Upon ligand binding, EGFR dimerizes, autophosphorylates and becomes active. Many signaling pathways are initiated through the EGFR and related ligands. Phosphorylation of Tyr 992 results in the binding and activation of Phospholipase-C gamma (PLC-gamma).

**Immunogen:** This EGF Receptor antibody was raised against a synthetic phosphopeptide corresponding to the amino acid residues region that contains Tyr992 of human Phosphorylated-EGFR.

**Buffer:** 12 mM sodium borate, 50 mM boric acid, 37 mM NaCl, 5 mM EDTA and 50% glycerol.

### Application Notes:

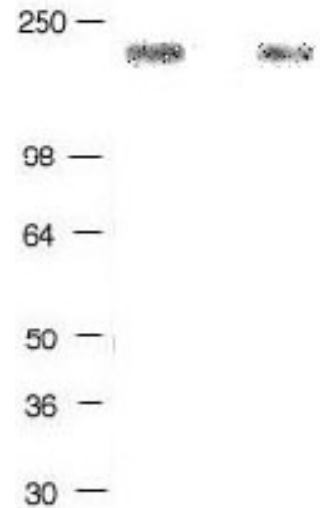
Applications Validated by Active Motif:

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



### EGF Receptor phospho Tyr992 pAb tested by Western blot.

Detection of Phospho-EGFR by Western blot. The analysis was performed using A-431 nuclear extract (EGF treated) and EGF Receptor phospho Tyr992 pAb at a 1:500 dilution. The reactions were incubated in the absence (lane 1) and presence of immunizing peptide (lane 2), and in the presence of non-phosphorylated peptide (lane 3), which confirms the specificity of the antibody for the phosphorylated form of the EGF receptor.