

## Recombinant ErbB-4 protein

**Catalog No:** 31167

**Lot No:** 02606001

**Expressed In:** Baculovirus

**Quantity:** 10 µg

**Concentration:** 0.232 µg/µl

**Source:** Human

**Buffer Contents:** 10 µg of Recombinant ErbB-4 protein in 50 mM Tris-HCl, pH 8.0; 100 mM NaCl, 5 mM DTT, 15 mM reduced glutathione, 20% glycerol.

**Protein Details:** Recombinant human ErbB-2 protein was produced using baculovirus infected Sf9 cells. The protein was made against amino acids R676-V1307, accession number NM\_005235 and N-terminally fused to a GST-HIS6-Thrombin cleavage site. Purified by GSH-agarose affinity purification.

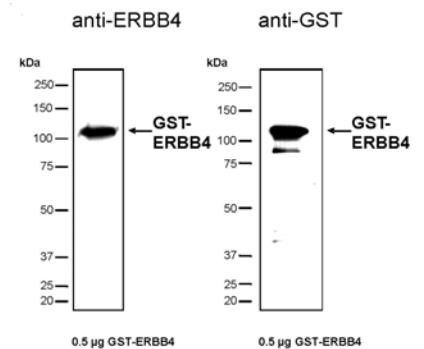
**Application Notes:** Recombinant ErbB-4 is suitable for kinase assays and Western blot. The molecular weight of the protein is ~102.062 kDa. The activity of the protein is ~ 26 pmol/µg min.

Recommended kinase reaction conditions: 60 mM HEPES-NaOH, pH 7.5, 3 mM MgCl<sub>2</sub>, 3 mM MnCl<sub>2</sub>, 3 µM Na-orthovanadate, 1.2 mM DTT, ATP (variable), 2.5 µg/50 µl PEG20.000, Substrate: PolyAEKY 5 µg/50 µl, Recombinant ErbB-4: 50 ng/50 µl.

Kinase activity may vary depending on the substrate and reaction conditions used.

**Storage and Guarantee:** Recombinant proteins in solution are temperature sensitive and must be stored at -80°C to prevent degradation. Avoid repeated freeze/thaw cycles and keep on ice when not in storage. This product is guaranteed for 6 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



### Recombinant ErbB-4 protein tested by SDS-PAGE gel.

0.5 µg of ErbB-4 protein was analyzed using SDS/PAGE followed by Western blot with either an anti-ErbB-4 or anti-GST antibody.