

## Anti-phospho-EPHB2 (Tyr469)

Rabbit Polyclonal Antibody

Catalog # E22-65R

Lot # J1274-14

### Cited Applications

WB

Suggested Dilutions:

WB 1:1,000

Ideal working dilutions for each application should be empirically determined by the investigator.

### Specificity

Recognizes the EPHB2 protein phosphorylated at tyrosine 469

### Cross Reactivity

Human, Mouse, Bovine, Canine, Chicken, non-Human Primate, Xenopus and Zebrafish

### Host/Isotype/Clone#

Rabbit, IgG

### Immunogen

Synthetic phospho-peptide corresponding to amino acid residues surrounding Tyr469 conjugated to KLH

### Formulation

100 µl in 10 mM HEPES (pH 7.5), 150 mM NaCl, 100 µg per ml BSA and 50% glycerol.

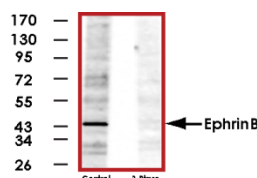
### Scientific Background

EPHB2 is a member of the Eph family of receptor tyrosine kinases that mediate neuro-developmental processes such as boundary formation, axon guidance, vasculogenesis and cell migration. Activated EPHB2 regulates the activity of the non-receptor tyrosine kinase Abl (1). EPHB2 is overexpressed in a number of tumors, particularly glioblastoma. In fact, EPHB2 overexpression leads to increased glioma cell migration and invasion (2).

### References

1. Yu, H H. et al: Multiple signaling interaction of Abl and Arg kinases with the EphB2 receptor. *Oncogene*. 2001 Jul 5;20(30):3995-4006.
2. Nakada, M. et al: The phosphorylation of EphB2 receptor regulates migration and invasion of human glioma cells. *Cancer Res*. 2004 May 1;64(9):3179-85.

### Sample Data



Western blot of rat testes lysate showing specific immunolabeling of the ~46kDa EPHB2 phosphorylated at Tyr469 (Control). The phosphospecificity of this labeling is shown in the second lane (lambda-phosphatase: λ-Ptase). The blot is identical to the control except that it was incubated in λ-Ptase (1200 units for 30 min) before being exposed to the Anti-phospho-EPHB2 (Tyr469). The immunolabeling of the EPHB2 band is completely eliminated by treatment with λ-Ptase.

## Anti-phospho-EPHB2 (Tyr469)

Rabbit Polyclonal Antibody

Catalog Number

E22-65R

Specific Lot Number

J1274-14

Purification Affinity chromatography

Stability 1yr at -20°C from date of shipment

Storage & Shipping

Store product at -20°C. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on ice packs.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: [orders@signalchem.com](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

**FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.**