

## Anti-phospho-ERK1/2 (Tyr204)

Rabbit Polyclonal Antibody

**Catalog # M29-65BR**

Lot # J1178-28

### Cited Applications

WB, IHC, IF

*Suggested Dilutions:*

WB: 1:500-1:1000

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

### Specificity

Recognizes the ERK1/2 protein phosphorylated at tyrosine 204

### Cross Reactivity

Human, Mouse and Rat

### Host/Isotype/Clone#

Rabbit, IgG

### Immunogen

Synthetic phospho-peptide corresponding to amino acid residues surrounding Tyr204

### Formulation

PBS (pH 7.4) 150mM NaCl, 0.02% sodium azide and 50% glycerol.

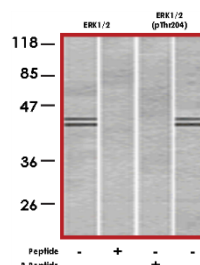
### Scientific Background

Extracellular-Signal Regulated Kinase/Mitogen-Activated Protein Kinase (ERK/MAPK) is a serine threonine kinase. It plays an integral role of cellular signaling during mitogenesis and differentiation of mitotic cells. ERK is presumed to have a key role in learning and memory (1,2,3). The activity of this kinase is regulated by phosphorylation at Thr202 and Tyr204 (4). Activated ERK1/2 translocates into the nucleus where it phosphorylates various transcription factors (e.g Elk-1, c-Myc, c-Jun, c-Fos and C/EBP beta).

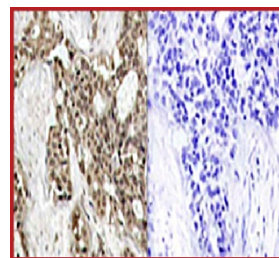
### References

1. Adams, J P. et al: Molecular psychology: Roles for the ERK MAP kinase cascade in memory. *Annu Rev Pharmacol Toxicol* 2002 42:135-163.
2. Johnson, G L. et al: Mitogen-activated protein kinase pathways mediated by ERK, JNK, and p38 protein kinases. *Science* 202 298:1911-1912.
3. Tanoue, T J. et al: Molecular recognitions in the MAP kinase cascades. *Cellular Signaling* 2003 15:455-462.
4. Ahn, N G.: The MAP kinase cascade. Discovery of a new signal transduction pathway. *Mol Cell Biochem* 1993 127-128:201-209.

### Sample Data



Western blots from NIH-3T3 cells using ERK1/2 antibody (lanes 1 and 2) and anti-phospho-ERK1/2 (Tyr204) antibody (lanes 3 and 4).



P-Peptide - +

Immunohistochemical analysis of paraffin-embedded human breast carcinoma tissue using anti-phospho-ERK1/2 (Tyr204) antibody.

## Anti-phospho-ERK1/2 (Tyr204)

Rabbit Polyclonal Antibody

Catalog Number

M29-65BR

Specific Lot Number

J1178-28

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.**