

## Anti-phospho-TRKA (Ser791)

Rabbit Polyclonal Antibody

**Catalog # N16-65R**

Lot # Z2014-43

### Cited Applications

IF

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

### Specificity

Recognizes the TRKA protein phosphorylated at serine 791

### Cross Reactivity

Human

### Host/Isotype/Clone#

Rabbit, IgG

### Immunogen

The antibody was produced against synthesized phosphopeptide derived from human TrKA around the phosphorylation site of tyrosine791 (P-V-YP-L-D).

### Formulation

Phosphate buffered saline (without Mg<sup>2+</sup> and Ca<sup>2+</sup>), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.

### Stability

1yr at -20°C from date of shipment

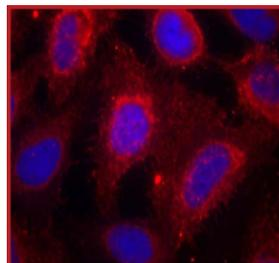
### Scientific Background

TRKA is a member of the TRK proto-oncogene family and encodes a 140-kilodalton, membrane-spanning protein tyrosine kinase that is the functional receptor for nerve growth factor (NGF). NGF elicits the rapid phosphorylation of gp140trk on tyrosine residues leading to increased c-Fos expression, DNA synthesis and morphologic transformation (1). A decreased expression of TRKA on the striatal cholinergic neurons has been observed which may contribute, when it reaches a crucial threshold, to the death of cholinergic neurons observed in Alzheimer disease (2).

### References

1. Kaplan, D R. et al: The trk proto-oncogene product: a signal transducing receptor for nerve growth factor. Science. 1991 Apr 26;252(5005):554-8.
2. Boissiere, F. et al: Neurotrophin receptors and selective loss of cholinergic neurons in Alzheimer disease. Mol Chem Neuropathol. 1996 May-Aug;28(1-3):219-23.

### Sample Data



Immunofluorescence staining of methanol-fixed HeLa cells using Anti-phospho-TRKA (Ser791).

## Anti-phospho-TRKA (Ser791)

Rabbit Polyclonal Antibody

Catalog # N16-65R

Lot # Z2014-43

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.**