

## Anti-phospho-EIF2AK2 (Thr446)

Rabbit Polyclonal Antibody

**Catalog # P80-65R**

Lot # J1178-14

### Cited Applications

WB, IHC

*Suggested Dilutions:*

WB: 1:500-1:1000 IHC: 1:50-1:100

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

### Specificity

Recognizes the EIF2AK2 protein phosphorylated at threonine 446

### Cross Reactivity

Human and Mouse

### Host/Isotype/Clone#

Rabbit, IgG

### Immunogen

Synthetic phospho-peptide corresponding to amino acid residues surrounding Thr446

### Formulation

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

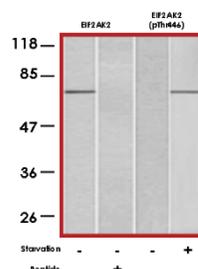
### Scientific Background

EIF2AK2 (also known as double-stranded RNA-activated protein kinase and PKR) is a protein kinase that is involved in HIV/gp120-associated neurodegeneration (1). EIF2AK2 is a mediator of gp120 neurotoxicity and is a substrate for a family of protein kinases involved in environmental stress responses. Activation of EIF2AK2 by dsRNA leads to autophosphorylation at Thr446/451. Activated EIF2AK2 phosphorylates the alpha subunit of EIF2, which in turn inhibits protein synthesis. EIF2AK2 plays a critical role in mRNA translation, cell proliferation and apoptosis. Crosstalk between EIF2AKs and p53 has implications in cell proliferation and tumorigenesis (2).

### References

1. Baltzis, D. et al: The eIF2alpha kinases PERK and PKR activate glycogen synthase kinase 3 to promote the proteasomal degradation of p53. *J. Biol Chem.* 2007; 282(43):31675-87
2. Alirezai, M. et al: Human immunodeficiency virus-1/surface glycoprotein 120 induces apoptosis through RNA-activated protein kinase signaling in neurons. *J. Neurosci.* 2007;27(41):11047-55.

### Sample Data



Western blot analysis of extracts from K562 cells using EIF2AK2 antibody (lanes 1 and 2) and anti-phospho-EIF2AK2 (Thr446) antibody (lanes 3 and 4).

## Anti-phospho-EIF2AK2 (Thr446)

Rabbit Polyclonal Antibody

Catalog Number

P80-65R

Specific Lot Number

J1178-14

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.**