

Anti-EIF2AK2 (PKR)

Rabbit Polyclonal Antibody

Catalog # **P80-63R**

Lot # M168-2

Cited Applications

- Western blot (1:1000)

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

Specificity

Recognizes the EIF2AK2 (PKR) protein

Cross Reactivity

- Western blot human EIF2AK2

EIF2AK2 from other species may also be detectable

Host

Rabbit

Immunogen

Synthetic peptide corresponding to amino acids 449-453 of human EIF2AK2

Formulation

PBS, pH7.4, 0.02%NaN₃, 50% glycerol

Stability

Store at 4°C (add 0.1% NaN₃) for several months, and at -20°C for longer periods. 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.

Scientific Background

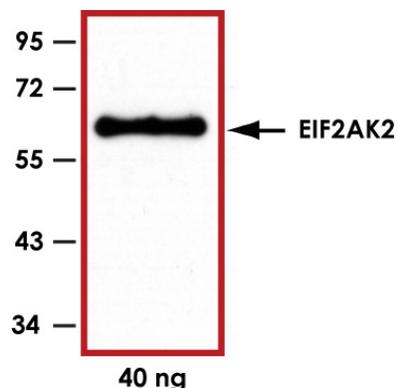
EIF2AK2 (also known as double-stranded RNA-activated protein kinase) has been shown to be involved in HIV/gp120-associated neurodegeneration (1). EIF2AK2 acts as a critical mediator of gp120 neurotoxicity and is a substrate for a family of protein kinases that respond to various forms of environmental stress. Activation of EIF2AK2 leads to its autophosphorylation and then phosphorylation of its natural substrate, the alpha subunit of eukaryotic protein synthesis initiation factor-2. A novel cross-talk between the EIF2AKs and p53 has been shown that has implications in cell proliferation and tumorigenesis (2).

References

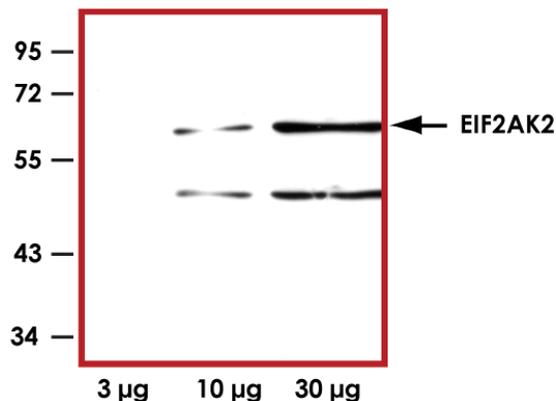
1. Baltzis, D. et al: The eIF2alpha kinases PERK and PKR activate glycogen synthase kinase 3 to promote the proteasomal degradation of p53. *J. Biol Chem.* 2007; 282(43):31675-87.
2. Alirezaei, M. et al: Human immunodeficiency virus-1/surface glycoprotein 120 induces apoptosis through RNA-activated protein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

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

Sample Data



Representative Western blot with Anti-EIF2AK2 (1:1000) using 40ng of GST-tagged human recombinant EIF2AK2.



Representative Western blot with Anti-EIF2AK2 (1:1000) using 3 µg, 10 µg and 30 µg of Jurkat cell lysate.

Anti-EIF2AK2 (PKR)

Rabbit Polyclonal Antibody

Catalog #	P80-63R
Lot #	M168-2
Purification	Affinity chromatography
Concentration	1.0 µg/µL
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.