

Anti-eIF4E

Rabbit Polyclonal Antibody

Catalog # **E34-63R**

Lot # M190-5

Cited Applications

- Western blot (1:1000)

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

Specificity

Recognizes the eIF4E protein

Cross Reactivity

- Western blot eIF4E from human and mouse cells
eIF4E from other species may also be detectable

Host

Rabbit

Immunogen

Synthesized peptide corresponds to amino acid 209 of human eIF4E protein.

Formulation

PBS, PH7.4, 0.02% NaN₃, 50% glycerol

Stability

Store at 4°C 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

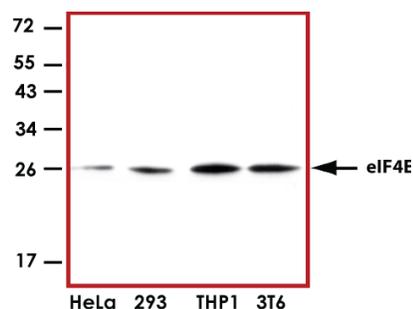
EIF4E is a eukaryotic translation initiation factor involved in directing ribosomes to the cap structure of mRNAs. EIF4E and is the rate-limiting component in protein synthesis. Overexpression of EIF4E causes malignant transformation (1). Overexpression of EIF4E in transgenic mice causes a marked increase in tumorigenesis. EIF4E is a potential physiological substrate for MNK1 and MNK2. MNK1 interacts with the EIF4F complex via its interaction with the C-terminal region of EIF4G, not through a direct interaction with EIF4E (2).

References

1. Ruggero D, et al: The translation factor eIF-4E promotes tumor formation and cooperates with c-Myc in lymphomagenesis. *Nature Med.* 10: 484-486, 2004.
2. Waskiewicz A J, et al: Mitogen-activated protein kinases activate the serine/threonine kinases Mnk1 and Mnk2. *EMBO J.* 16: 1909-1920, 1997.

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-eIF4E (1:1000) using 20 µg of HeLa, 293, THP1 and 3T6 cell lysates.

Anti-eIF4E

Rabbit Polyclonal Antibody

Catalog Number

E34-63R

Specific Lot Number

M190-5

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.