elF4E-BP2 polyclonal antibody

The eukaryotic translation initiation factors (eIFs) regulate translation at the level of messenger RNA engagement by the ribosome. eIF-2alpha promotes the binding of the initiator tRNA to 40S ribosomal subunits, and is a key signaling component of the unfolded protein response (UPR). eIF-4E recruits mRNAs to the eIF4F complex through binding of their 5 N7-methylguanosine cap. eIF-4E binding proteins-1 and -2 are translational repressors, sequestering eIF-4E in conditions of low cellular demand for translation. Their dissociation from eIF-4E is regulated by phosphorylation in response to growth-promoting stimuli.

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

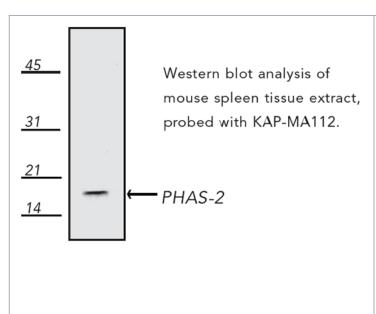
Order Online »

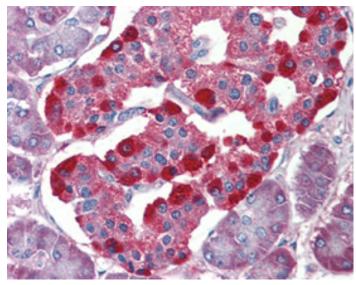
ADI-KAP-MA112-E

100µg

Manuals, SDS & CofA

View Online »





IHC Analysis: Human pancreas tissue stained with eIF4E-BP2, pAb at 10µg/ml.

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Eukaryotic translation initiation factor 4E-binding protein 2

Application IHC (PS), WB

Formulation Liquid. In borate buffered saline, pH 8.4, containing 25mM

sodium borate, 100mM boric acid, 75mM sodium chloride,

and 5mM EDTA.

GenBank ID L36056

Host Rabbit

Immunogen Synthetic peptide corresponding to a portion of human

eIF4E-BP2. The sequence is completely conserved in

mouse.

Purity Detail Peptide affinity purified.

Recommendation Dilutions/Conditions Western Blot (4µg/ml colorimetric, 4µg/ml ECL)Suggested

dilutions/conditions may not be available for all applications. Optimal conditions must be determined

individually for each application.

Source Purified from rabbit serum.

Species Reactivity Bovine, Dog, Human, Mouse, Porcine, Rabbit, Rat

UniProt ID Q13542

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

•

Last modified: May 29, 2024



info-

eu@enzolifesciences.com