

eIF4E-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

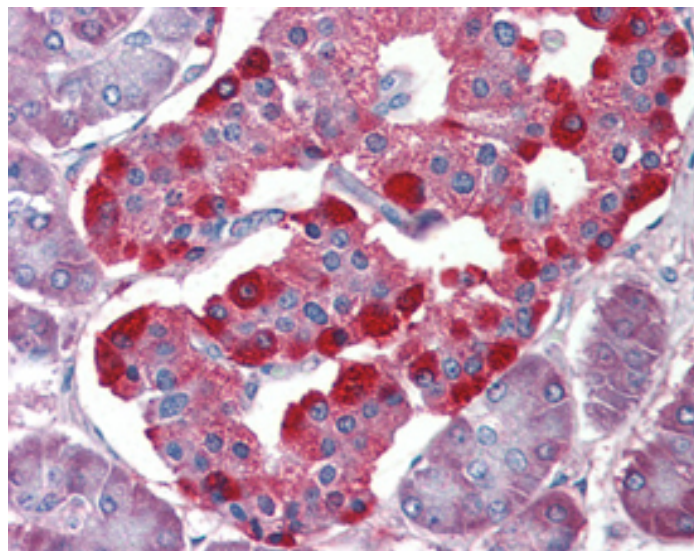
[View Online »](#)

45
31
21
14



Western blot analysis of
mouse spleen tissue extract,
probed with KAP-MA112.

← *PHAS-2*



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 .



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)