

eIF2 α (human), (recombinant) (His- tag)

Recombinant human eIF2 α (His-tag) that regulates translation initiation and stress-

induced signaling.
Eukaryotic translation initiation factor 2 α (eIF2 α) promotes the binding of initiator tRNA to the 40S ribosomal subunits, and controls translational rates via phosphorylation/dephosphorylation mechanisms. The double-stranded RNA (dsRNA)-activated protein kinase (PKR) provides an important control in the regulation of protein synthesis initiation through the phosphorylation of eIF2 α . Phosphorylated eIF2 α inhibits cell growth and induces apoptosis. Phosphorylation of eIF2 α occurs under a variety of conditions, including viral infection, apoptosis, nutrient deprivation, heme-deprivation, and certain stresses. eIF2 exists as a heterotrimeric protein composed of 3 nonidentical subunits: α (36kDa), β (38kDa), and γ (52kDa).

Citations: 3

[View Online »](#)

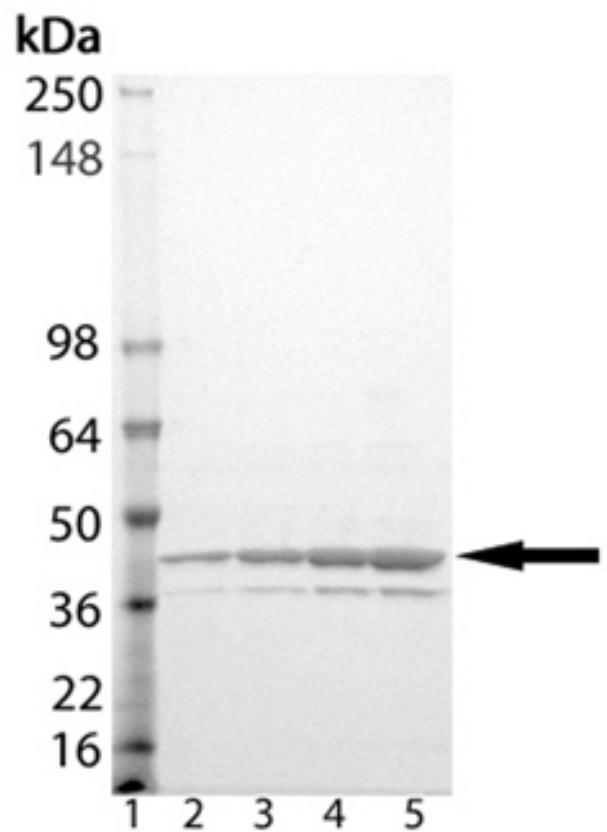
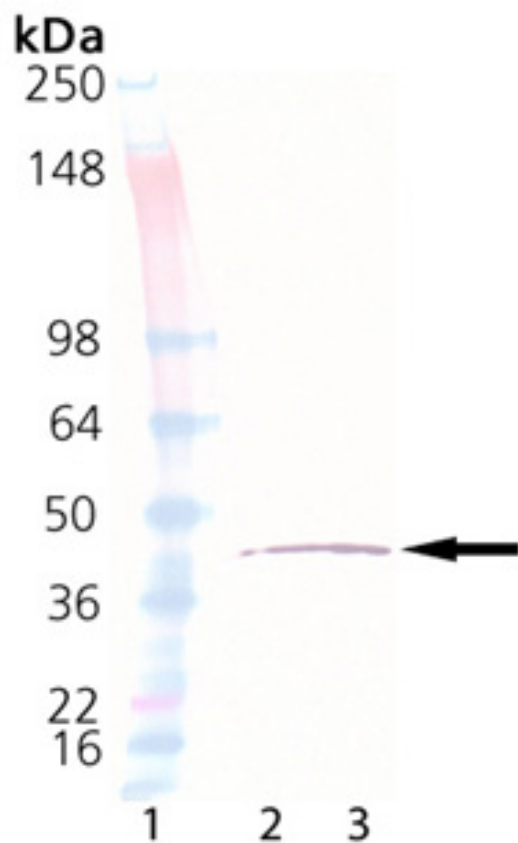
Ordering Information

[Order Online »](#)

ADI-KPR-CP132-0050	50 μ g
ADI-KPR-CP132-0200	200 μ g

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name Eukaryotic translation initiation factor 2 subunit α

Application Notes Western blot control.

Formulation Liquid. In 50mM TRIS, pH 7.5, containing 10% glycerol, 0.5% Triton X-100, 300mM potassium chloride, and 4mM DTT.

MW ~36kDa

Purity $\geq 85\%$ (SDS-PAGE; Western blot)

Purity Detail Purified by multi-step chromatography.

Source Produced in *E. coli*. Human eIF2 α is fused at the N-terminus to a His-tag.

UniProt ID P05198

Last modified: October 9, 2025



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)