

[pSer⁵²]Eukaryotic Initiation factor 2 α polyclonal antibody

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 6

[View Online »](#)

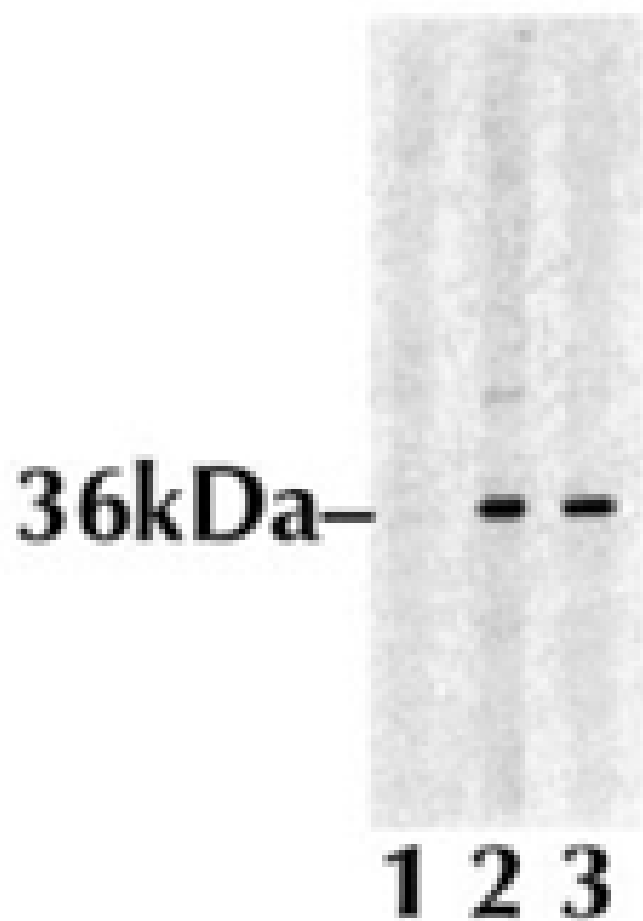
Ordering Information

[Order Online »](#)

BML-SA405-0020	20 μ l
BML-SA405-0100	100 μ l

Manuals, SDS & CofA

[View Online »](#)



Mouse 3T3 cell extracts were blotted and incubated with anti-eIF2- α (pS52) antibody following (1) prior incubation in the presence of the peptide immunogen, (2) the nonphosphopeptide corresponding to the eIF2- α (pS51) phosphopeptide, (3) or no peptide. Only the phospho-peptide corresponding to the region surrounding serine 51 (serine 52 in mouse) blocks the antibody signal.

Handling & Storage

Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -80°C.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	eIF2α
Application	WB
Crossreactivity	Does not cross-react with the unphosphorylated form of eIF2α.
Formulation	Liquid. In Dulbecco's PBS, pH 7.3, containing 1mg/ml BSA, 50% glycerol and 0.05% sodium azide.
Host	Rabbit
Immunogen	Synthetic peptide corresponding to a portion of human eIF2α (eukaryotic initiation factor-2α) phosphorylated at Ser51(Ser52 in mouse).
Recommendation Dilutions/Conditions	Western Blot (1:1,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human, Mouse, Yeast
Specificity	Recognizes human and yeast eIF2α phosphorylated at Ser ⁵¹ and mouse eIF2α phosphorylated at Ser ⁵² .
UniProt ID	P05198

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-be@enzolifesciences.com)

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

Last modified: May 29, 2024

For local distributors and detailed product information visit us online: enzo.com All product names, logos, and brands are the property of their respective owners.