

Actin clonal antibody (S12-I)

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

Citations: 1

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

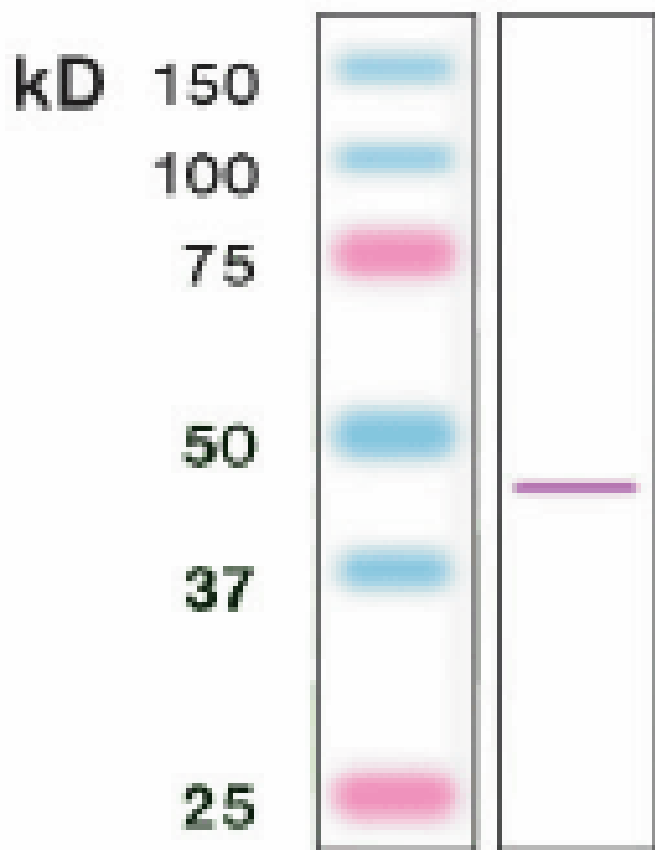


Figure: Western blot of actin in mouse brain crude lysate (50 μ g of protein loaded).

Handling & Storage

Use/Stability	Stable for at least 12 months after receipt when stored as recommended.
Handling	Avoid freeze/thaw cycles.
Long Term Storage	+4°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application	ELISA, IP, WB
Clone	S12-I
Formulation	Liquid. In 20mM TRIS/HCl, pH 8.0, containing 20mg/ml BSA and 0.05% sodium azide.
Host	Rabbit
Immunogen	Synthetic peptide corresponding to a C-terminal sequence of human β -actin.
Isotype	IgG
Purity Detail	Major clone of IgGs obtained from immunoaffinity purified immunoglobulins corresponding to immunogenic peptide.
Recommendation Dilutions/Conditions	ELISA (1:10,000-1:20,000)Western Blot (1:2,000; Wash buffer: 1x TRIS buffered saline (TBS) containing 0.1% Triton X-100, Blocking buffer: 1x TBS containing 0.1% Triton X-100 and 5% nonfat dry milk. Incubate the membrane with antibody diluted in blocking buffer for 2 hours at room temperature)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human, Mouse, Rat
Specificity	Recognizes α -actin and β -actin.

Clonal Rabbit Antibody is a pure homogenous fraction of rabbit immunoglobulin (IgG) which corresponds to a unique linear epitope on the target protein. This epitope is designed after a detailed analysis of the antigen structure, regulatory post-translational modifications, protein/protein interaction domain and protein folding. The chosen sequence (peptide) is synthesized and the rabbit is immunized. The clone is then separated from the crude antiserum by “*in vitro* cloning technology”. The resulting product is monospecific and characterized by superior specificity, affinity and avidity.

UniProt ID

P60709 (β-actin)

Worry-free Guarantee

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

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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