

Leptin receptor polyclonal antibody

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

Citations: 1

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

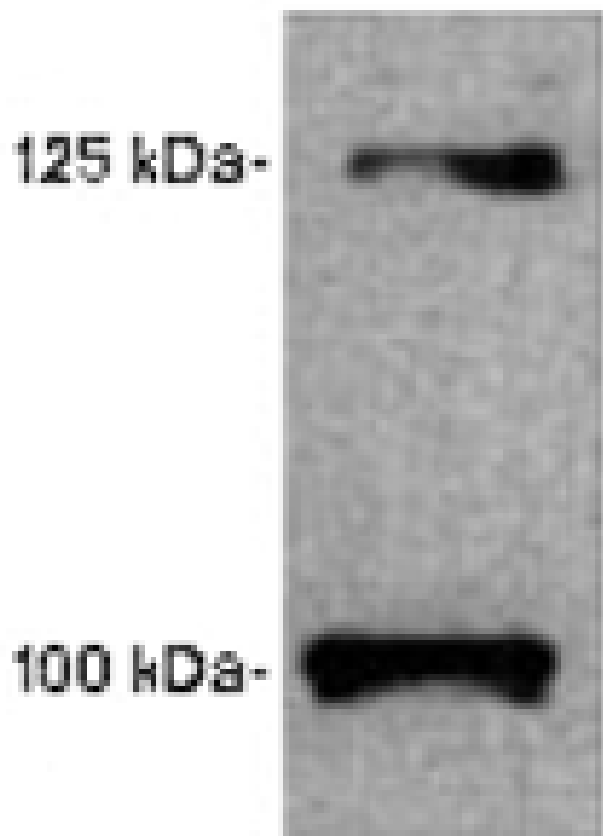


Figure: Western blot of both short and long forms of leptin receptor on mouse lung tissue extract using PAb to Leptin Receptor ([ALX-210-168](#)).¹, Leptin receptor polyclonal antibody Western blot, Leptin receptor polyclonal antibody Western blot, yes ALX-210-169, 210-169.gif, <http://static.enzolifesciences.com/fileadmin/files/image/210-169.gif>, **Figure:** Immunolocalization of IP₃R-I in rat cerebellum using PAb to Inositol 1

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	OB-R, CD295
Application	WB
Application Notes	Detects bands of ~100kDa and ~125kDa by Western blot, representing the short and long forms of leptin receptor. Also detects two unrelated bands of ~31kDa and ~35kDa.
Formulation	Liquid. In PBS containing 0.05% sodium azide.
Host	Rabbit
Immunogen	Synthetic peptide corresponding to aa 577-594 (K577EIQWKTHEVFDKSKSA594) of the extracellular domain of rat leptin receptor.
Purity Detail	Purified IgG.
Recommendation Dilutions/Conditions	Western Blot (1:2,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human, Mouse, Rat
UniProt ID	Q62959

Worry-free Guarantee

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



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

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

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

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

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

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

Last modified: May 29, 2024