

# GLP1R polyclonal antibody

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

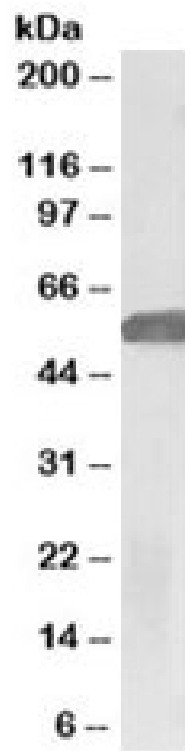
## Ordering Information

[Order Online »](#)

ENZ-ABS892-0100	100µg
-----------------	-------

## Manuals, SDS & CofA

[View Online »](#)



Western blot testing of human hippocampus lysate with GLP1R antibody at 1 $\mu$ g/ml. Predicted molecular weight ~53kDa.

## Handling & Storage

**Handling** Aliquot and store the GLP1R antibody at -20°C.

**Long Term Storage** -20°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** Glucagon-like peptide 1 receptor

**Application** ELISA, WB

**Formulation** Liquid. 0.5 mg/ml in 1X TBS, pH7.3, containing 0.5% BSA (US sourced) and 0.02% sodium azide.

**Host** Goat

**Immunogen** Amino acid sequence from the extracellular portion of GL1PR (QHQWDGLLSYQD).

**Isotype** Ig

**Purity Detail** Antigen affinity purified.

**Recommendation Dilutions/Conditions** ELISA (peptide) (LOD: 1:128,000)Western blot (1-3µg/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

**Species Reactivity** Human

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

Last modified: July 12, 2024



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-  
fr@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)