

# GLP-1 (7-37, 7-36, amide, free NT) monoclonal antibody (4)

Glucagon-like peptide 1 (7-36) amide (GLP-1 (7-36) amide) is the principal active form of GLP-1, the other being GLP-1 (7-37). GLP-1 is a peptide hormone of the glucagon family, produced by the L cells of the intestinal mucosa from the same prohormone as glucagon. The active forms are potent stimulators of glucose-dependent insulin secretion. The sequence of GLP-1 is fully conserved in all mammalian species examined so far.

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

Citations: 1

[View Online »](#)

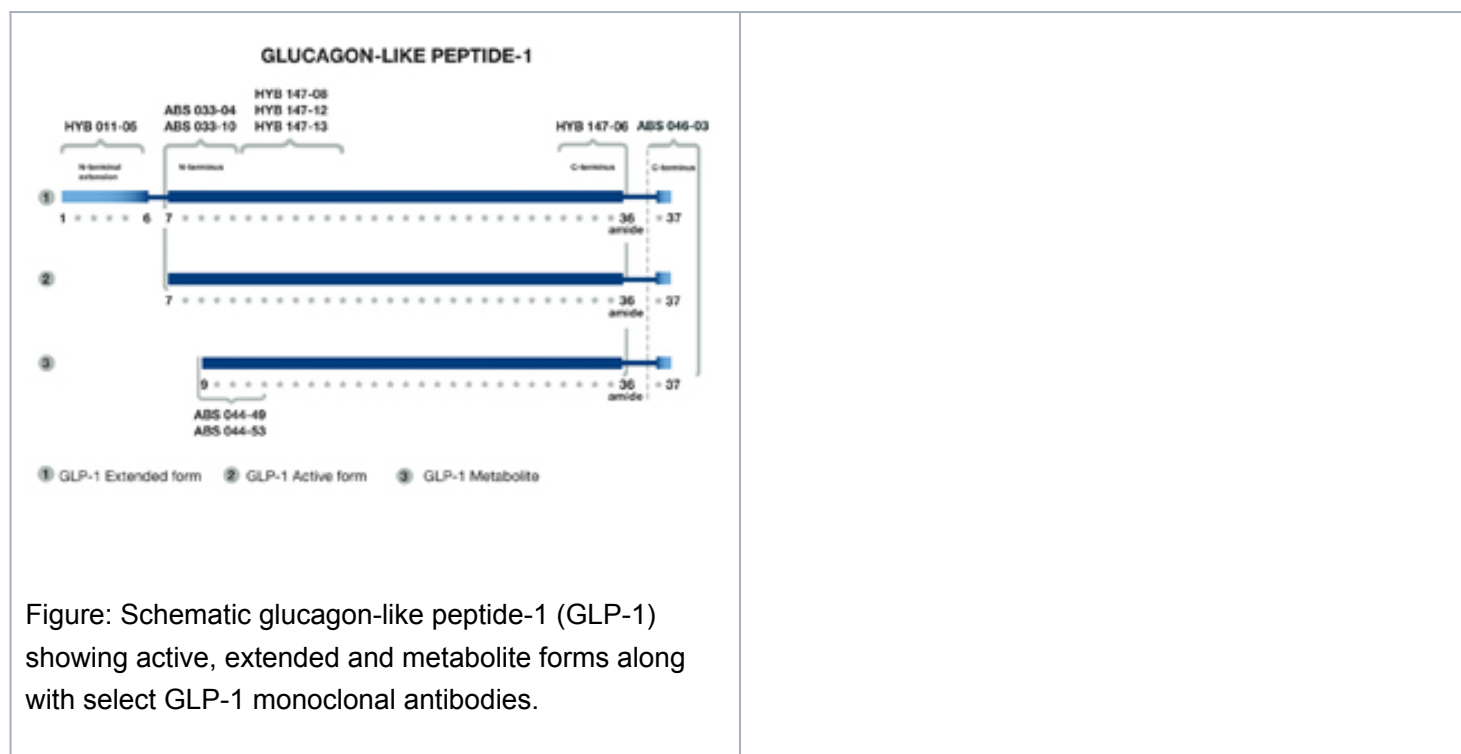
## Ordering Information

[Order Online »](#)

BPD-ABS-033-04-1	1mg
BPD-ABS-033-04-04	400µg

## Manuals, SDS & CofA

[View Online »](#)



# Handling & Storage

Short Term Storage +4°C

Shipping Blue Ice

# Regulatory Status

RUO - Research Use Only

# Product Details

Alternative Name	Glucagon like peptide 1
Application	ELISA, IHC
Application Notes	Not recommended for use in sandwich ELISA.
Clone	4
Crossreactivity	Shows <0.2% cross-reactivity with GLP-1(1-37), GLP-1(9-36)amide, glucagon, human GIP and exendin-4. Cross-reacts approximately 1% with human GLP-2.
Formulation	Liquid. In 0.01M phosphate buffer, pH 7.4, containing 0.5M sodium chloride and 15mM sodium azide.
Host	Mouse
Immunogen	Synthetic peptide corresponding to aa 7-17 of GLP-1.
Isotype	IgG1κ
Purity Detail	Protein A/G-affinity purified.
Recommendation Dilutions/Conditions	Immunohistochemistry (1:500)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Mammalian
UniProt ID	P01275
Worry-free Guarantee	This antibody is covered by our <a href="#">Worry-Free Guarantee</a> .



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)