

# GLP-1 (7-37, 7-36, amide, free NT) monoclonal antibody (10) (biotin conjugate)

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](#).

## Ordering Information

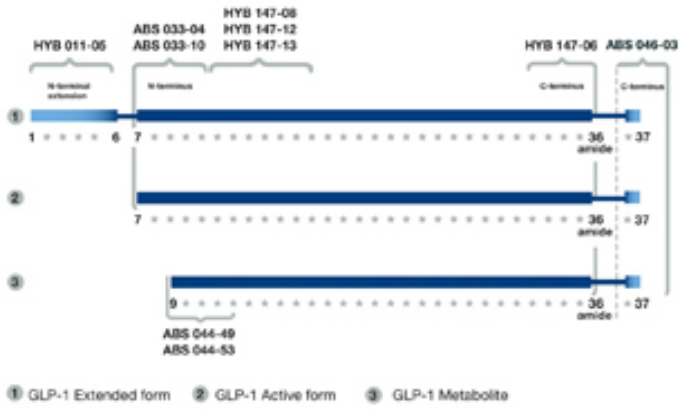
[Order Online »](#)

BPD-ABS-033-10B-015	150µg
---------------------	-------

## Manuals, SDS & CofA

[View Online »](#)

# GLUCAGON-LIKE PEPTIDE-1



Schematic glucagon-like peptide-1 (GLP-1) showing active, extended and metabolite forms along with select GLP-1 monoclonal antibodies.

## 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

Clone 10

**Crossreactivity** Shows <0.2% cross-reactivity with GLP-1(9-36)amide, glucagon and human GIP. Cross-reacts 0.25% with GLP-1(1-37), approximately 6% with human GLP-2 and approximately 50% with exendin-4.

**Formulation** Liquid. In 0.01M phosphate buffer, pH 7.4, containing 0.14mM sodium chloride and 15mM sodium azide.

Host Mouse

**Immunogen** Synthetic peptide corresponding to aa 7-17 of GLP-1.

Isotype IgG2ak

**Purity Detail** Protein A/G-affinity purified.

**Species Reactivity** Mammalian

UniProt ID P01275

**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-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)