

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

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

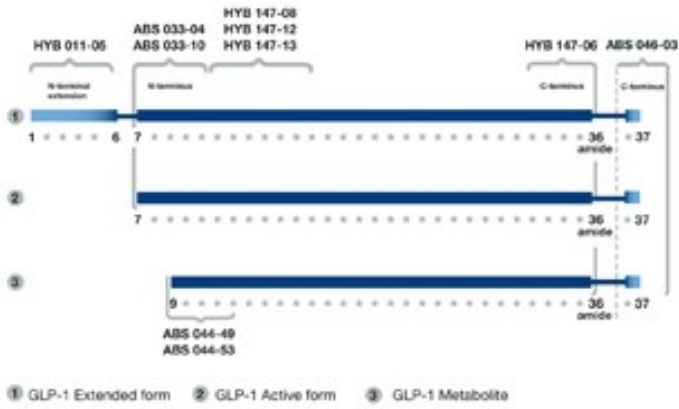
[Order Online »](#)

BPD-ABS-033-10-1	1mg
BPD-ABS-033-10-04	400µg

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

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Glucagon like peptide 1
Application	ELISA, WB
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 0.5M 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