GLP-1 (9-37, 9-36, amide, free NT) monoclonal antibody (49)

Glucagon-like peptide-1 (9-36) amide and GLP-1 (9-37) are the forms of GLP-1 which result from rapid degradation of the active forms of the peptide (GLP-1 (7-36) amide and GLP-1 (7-37)) by the enzyme dipeptidyl peptidase-IV (DPP-IV, also known as CD26 or adenosine deaminase binding protein). 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-044-49-1	1mg
BPD-ABS-044-49-04	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, WB

Clone 49

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 9-18 of GLP-1.

lgG1κ

Purity Detail Protein A-affinity purified.

Species Reactivity Mammalian

UniProt ID P01275

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

.

Last modified: May 29, 2024

