GLP-1 (Arg-Gly, nonamidated, CT) monoclonal antibody (03)

Glucagon-like peptide-1 (GLP-1) is a peptide hormone of the glucagon family released in the intestine in response to eating. GLP-1 is responsible for maintaining the correct balance between blood glucose levels and levels of glucose being adsorbed in the intestine. The active forms of GLP-1 are potent stimulators of glucose-dependent insulin secretion and may be drug candidates for the treatment of obesity and type 2 diabetes.

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

Order Online »

BPD-ABS-046-03-1	1mg
BPD-ABS-046-03-04	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 03

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 9-37 of GLP-1.

lgG1κ

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

.

Last modified: May 29, 2024

