

GLP-1 (Arg-Gly, non-amidated, 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-046-03-04	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.
Isotype	IgG1κ
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



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)