

Glucagon monoclonal antibody (09)

Glucagon is a peptide hormone of 29 amino-acid residues produced by the alpha cells of the pancreatic islets. Its effect is to raise blood sugar as a stress hormone and counter-regulatory hormone for insulin. It is produced by cell-type-specific processing of proglucagon, which in the alpha cells is split to form glicentin-related pancreatic polypeptide, glucagon and major proglucagon fragment, while the same precursor is split to form glicentin, oxyntomodulin (C-terminally extended glucagon-37), GLP-1 and GLP-2 in the L cells of the intestinal mucosa.

This antibody is covered by our [Worry-Free Guarantee](#).

Ordering Information

[Order Online »](#)

BPD-ABS-061-09-1	1mg
BPD-ABS-061-09-04	400µg

Manuals, SDS & CofA

[View Online »](#)

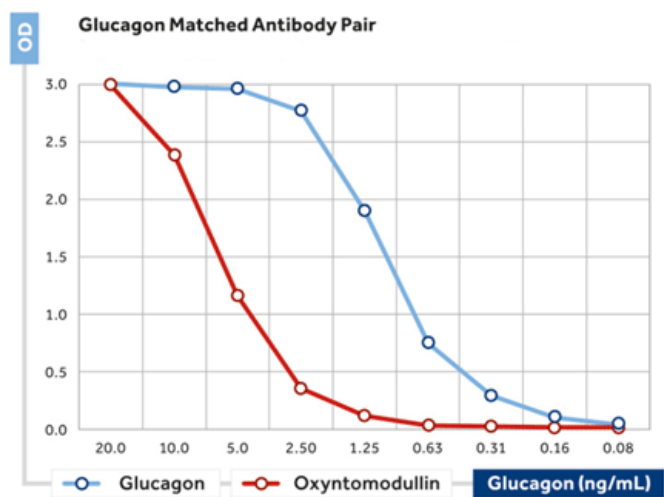


Figure 1: ELISA. Calibration curve for matched antibody pair Glucagon, mAb (09) (Prod. No. BPD-ABS-061-09) and Glucagon, mAb (08) (biotin conjugate) (Prod. No. BPD-ABS-058-08B)

Handling & Storage

Short Term Storage +4°C

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application ELISA

Clone 09

Formulation Liquid. In 0.01M phosphate buffer, pH 7.4, containing 0.5M sodium chloride and 15mM sodium azide.

Host Mouse

Immunogen Synthetic C-terminal fragment corresponding to human glucagon coupled to carrier.

Isotype IgG1κ

Purity Detail Protein A-affinity purified.

Species Reactivity Human, Porcine, Rat

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)