

# Glucagon monoclonal antibody (08) (biotin conjugate)

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 »](#)

<b>BPD-ABS-058-08B-015</b>	150µg
----------------------------	-------

## Manuals, SDS & CofA

[View Online »](#)

## 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	08
Formulation	Liquid. In 0.01M phosphate buffer, pH 7.4, containing 0.14M sodium chloride and 15mM sodium azide.
Host	Mouse
Immunogen	Synthetic N-terminal fragment corresponding to human glucagon coupled to carrier.
Isotype	IgG1k
Purity Detail	Protein A-affinity purified.
Species Reactivity	Human, Porcine, Rat
UniProt ID	P01275
Worry-free Guarantee	This antibody is covered by our <a href="#">Worry-Free Guarantee</a>

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-  
be@enzolifesciences.com](mailto:info-be@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)