

Oxyntomodulin monoclonal antibody (46)

Oxyntomodulin is a peptide hormone of 37 amino-acid residues (29-amino acid sequence of glucagon, with an 8-amino acid carboxy-terminal extension) produced by the L cells of the gut. It modulates gastric acid secretion from the oxyntic cells of the stomach, inhibits appetite and has glucose lowering effects. Oxyntomodulin is produced by cell-type-specific processing of proglucagon that forms as well glicentin, 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-064-46-1	1mg
BPD-ABS-064-46-04	400µg

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application ELISA

Clone 46

Crossreactivity Does not cross-react with glucagon.

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

GenBank ID 2641

Host Mouse

Immunogen Synthetic oxyntomodulin 31-37 coupled to carrier.

Isotype IgG1κ

Purity Detail Protein A/G-purified.

Source From culture supernatant.

Species Reactivity Human, Mouse, Rat

Specificity Specific for C-terminus of human oxyntomodulin.

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