

Oxyntomodulin (human) monoclonal antibody (20) (biotin conjugate)

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-059-20B-015	150µg
---------------------	-------

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 20

Crossreactivity Cross-reacts fully 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 full length oxyntomodulin coupled to carrier

Isotype IgG1κ

Purity Detail Protein A/G-purified.

Species Reactivity Human, Mouse, 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-
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)