

Oxyntomodulin monoclonal antibody (23)

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-064-23-1	1mg
BPD-ABS-064-23-04	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	23
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](#)